



Agilent SLIMS 6.7

Open Source Licenses

Agilent SLIMS contains software provided by third parties, including Open-Source software.
 This document contains licensing information about those third-party products used by Agilent SLIMS 6.7.

Title	Authors	Declared License
ActiveMQ :: Camel		Apache-2.0
ActiveMQ :: KahaDB Store		Apache-2.0
Apache Commons CLI	Beluga Behr, Peter Donald, Brian Egge, Duncan Jones, Berin Loritsch, Peter Maddocks, Alexandru Mocanu, Andrew Shirley, Greg Thomas, Slawek Zachcial, James Strachan, Bob McWhirter, John Keyes, Rob Oxspring, Emmanuel Bourg, Thomas Neidhart	Apache-2.0
Apache Commons Codec	Christopher O'Brien, Martin Redington, Jeffery Dever, Steve Zimmermann, Benjamin Walstrum, Oleg Kalnichevski, Dave Dribin, Alex Karasulu, Matthew Inger, Jochen Wiedmann, Sebastian Bazley, Matthew Pocock, Colm Rice, Henri Yandell, Tim OBrien, Scott Sanders, Rodney Waldhoff, Daniel Rall, Jon S. Stevens, Gary Gregory	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause

Apache Commons Codec	Christopher O'Brien, Martin Redington, Jeffery Dever, Steve Zimmermann, Benjamin Walstrum, Oleg Kalnichevski, Dave Dribin, Alex Karasulu, Matthew Inger, Jochen Wiedmann, Sebastian Bazley, Matthew Pocock, Colm Rice, Henri Yandell, Tim O'Brien, Scott Sanders, Rodney Waldhoff, Daniel Rall, Jon S. Stevens, Gary Gregory	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause
Apache Commons CSV	Henri Yandell, Martin van den Bernt, Yonik Seeley, Emmanuel Bourg, Gary Gregory, Benedikt Ritter, Rob Tompkins	Apache-2.0
Apache Commons CSV	Henri Yandell, Martin van den Bernt, Yonik Seeley, Emmanuel Bourg, Gary Gregory, Benedikt Ritter, Rob Tompkins	Apache-2.0
Apache Commons DBCP	Todd Carmichael, Wayne Woodfield, Dain Sundstrom, Philippe Mouawad, Glenn L. Nielsen, James House, James Ring, Peter Wicks, Morgan Delagrang, Geir Magnusson, Craig McClanahan, John McNally, Martin Poeschl, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Yoav Shapira, Jürg Schaible, Mark Thomas, Gary Gregory	Apache-2.0
Apache Commons Exec	Niklas Gustavsson, Benjamin Bentmann, Marco Ferrante, Jerome Lacoste, Milos Kleint, Pablo Hoertner, Niall Pemberton, Brett Porter, Trygve Laugstøl, Siegfried Goeschl, Sebastian Bazley, Gary Gregory	Apache-2.0

Apache Commons FileUpload	Aaron Freeman, Daniel Fabian, Gary Gregory, Jörg Heinicke, Stepan Koltsov, Michael Macaluso, Amichai Rothman, Alexander Sova, Paul Spurr, Thomas Vandahl, Henry Yandell, Jan Novotný, frank, Rafal Krzewski, Sean Legassick, Oleg Kalnichevski, David Sean Taylor, Martin Cooper, dlon Gillard, John McNally	Apache-2.0
Apache Commons IO	Rahul Akolkar, Jason Anderson, Nathan Beyer, Emmanuel Bourg, Chris Eldredge, Magnus Grimsell, Jim Harrington, Thomas Ledoux, Andy Lehane, Marcelo Liberato, Alban Peignier, Adam Retter, Ian Springer, Dominik Stadler, Masato Tezuka, James Urie, Frank W. Zammetti, Scott Sanders, dlon Gillard, Nicola Ken Barozzi	Apache-2.0
Apache Commons Logging		Apache-2.0
Apache Commons Text	Don Jeba, Sampanna Kahu, Jarek Strzelecki, Lee Adcock, Amey Jadiye, Arun Vinud S S, Ioannis Sermetziadis, Jostein Tveit, Luciano Medallia, Jan Martin Keil, Nandor Kollar, Nick Wong, Bruno P. Kinoshita, Benedikt Ritter, Rob Tompkins, Gary Gregory, Duncan Jones	Apache-2.0
Apache Commons Validator	Don Brown, Martin Cooper, David Graham, Ted Husted, Rob Leland, Craig McClanahan, James Mitchell, Niall Pemberton, James Turner, David Winterfeldt, Henri Yandell, Ben Speakmon, Nick Burch, Simone Tripodi, Benedikt Ritter, Gary Gregory	Apache-2.0
Apache FontBox		Apache-2.0; BSD-3-Clause
Apache Geronimo JMS Spec 2.0		Apache-2.0
Apache HttpClient Mime		Apache-2.0
Apache Log4j API		Apache-2.0; Multi-license: Apache-2.0 OR LGPL-3.0-only
Apache Log4j Core		Apache-2.0; Multi-license: Apache-2.0 OR LGPL-3.0-only
Apache Log4j SLF4J Binding		Apache-2.0
Apache Log4j Web		Apache-2.0

Apache PDFBox		Unicode; Zlib; BSD-3-Clause; APAFML; OFL-1.1; Apache-2.0
Apache PDFBox Debugger		Apache-2.0; Multi-license: AML OR Apache-2.0
Apache PDFBox tools		Apache-2.0
Apache POI		Apache-2.0; Multi-license: Apache-2.0 OR MIT
Apache POI		Apache-2.0
Apache POI		Apache-2.0
Apache Tika parser modules		Apache-2.0; Multi-license: Apache-2.0 OR LGPL-3.0-only
Apache Velocity	Will Glass-Husain, Geir Magnusson Jr., Daniel Rall, Henning P. Schmiedehausen, Nathan Bubna	Apache-2.0
ARQ		Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause; JSON; public-domain
ASM based accessors helper used by json-smart		Apache-2.0
AutoValue Annotations		Apache-2.0
barbecue	Sean C. Sullivan, Werner Keil	BSD-3-Clause
Barcode4J	Jeremias MÅrki, Web Maestro Clay Leeds, Nicola Ken Barozzi, Dietmar BÅ¼rkle	Apache-2.0
batik-awt-util		Apache-2.0
batik-bridge		Apache-2.0
batik-dom		Apache-2.0; W3C
batik-svggen		Apache-2.0
batik-util		Apache-2.0
batik-xml		Apache-2.0
Bean Validation API	Emmanuel Bernard, Hardy Ferentschik, Gunnar Morling	Apache-2.0

BeanUtils	Robert Burrell Donkin, dlon Gillard, Craig McClanahan, Geir Magnusson Jr., Scott Sanders, James Strachan, Rodney Waldhoff, Martin van den Bemt	Apache-2.0
Bouncy Castle Provider		Bouncy Castle Licence; public-domain; MIT
Bouncy Castle S/MIME API		Bouncy Castle Licence
Bouncy Castle TSP/MIME API		Bouncy Castle Licence
Camel :: Core		Apache-2.0
Camel :: HazelCast		Apache-2.0
Camel :: JMS		Apache-2.0
Camel :: Mail		Apache-2.0
Camel :: Netty		Apache-2.0
Camel :: Netty4		Apache-2.0
Camel :: Quartz2		Apache-2.0
Camel :: Spring		Apache-2.0
Camel :: Spring Integration		Apache-2.0
Camel :: Test		Apache-2.0
Camel :: Test :: Spring 4.1+		Apache-2.0
cglib		Multi-license: Apache-2.0 OR ASF 2.0; Apache-2.0; BSD-3-Clause
Commons DBCP	Todd Carmichael, Wayne Woodfield, Morgan Delagrange, Geir Magnusson, Craig McClanahan, John McNally, Martin Poeschl, Rodney Waldhoff, David Weinrich, Dirk Verbeeck	Apache-2.0
Commons JXPath	Dmitri Plotnikov, Craig McClanahan, Matt Benson	Apache-2.0

Commons Lang	Ringo De Smet, Steve Downey, Daniel Rall, Stephen Colebourne, Henri Yandell, Steven Caswell, Robert Burrell Donkin, Gary Gregory	Apache-2.0; Multi-license: Apache-2.0 OR GPL-2.0-only OR LGPL-2.1-only
Commons Math	C. Scott Ananian, Mikkel Meyer Andersen, Mark Anderson, RÅ@mi Arntzen, Michael Bjorkegren, John Bollinger, Cyril Briquet, Paul Cowan, Benjamin Croizet, Larry Diamond, Rodrigo di Lorenzo Lopes, Hasan Diwan, Ted Dunning, John Gant, Ken Geis, Bernhard GrÅ¼newaldt, Elliotte Rusty Harold, Matthias Hummel, Ismael Juma, Eugene Kirpichov	Apache-2.0; Multi-license: Apache-2.0 OR BSD-2-Clause; Multi-license: Apache-2.0 OR BSD-3-Clause
commons-httpclient	Armando Anton, Sebastian Bazley, Ola Berg, Sam Berlin, Mike Bowler, Samit Jain, Eric Johnson, Christian Kohlschuetter, Ryan Lubke, Sam Maloney, Rob Di Marco, Juergen Pill, Mohammad Rezaei, Roland Weber, Laura Werner, Mikael Wilstrom, Michael Becke, Jeff Dever, dlon Gillard, Ortwin Glueck	Apache-2.0; Multi-license: Apache-2.0 OR LGPL-2.0-or-later
Date Calc JODA		Apache-2.0
Date Calculator Common		Apache-2.0
Discovery	Richard Sitze, Craig R. McClanahan, Costin Manolache, James Strachan,	Apache-2.0

	Matthew Hawthorne, Davanum Srinivas, Rory Winston, Robert Burrell Donkin	
dom4j	Curt Arnold, David Lucas, David White, Jakob Jenkov, James Strachan, Laramie Crocker, Maarten Coene, Michael Skells, Nick Sanderson, Steen Lehmann, Tobias Rademacher, Bob McWhirter, Todd Wolff, OuYang Chen, Yuxin Ruan	BSD-3-Clause; SPL-1.0
EventBinder		Apache-2.0
FindBugs-jsr305		Apache-2.0; CC-BY-2.5
fr.opensagres.xdocreport.converter.docx.xwpf		MIT
GIN (GWT INjection)	Brian Stoler, Peter Schmitt, Jesse Wilson, Robbie Vanbrabant, Fazal Asim, Daniel Burrows, Scott Blum, Adam Winer, Philippe Beaudoin, Ben Chambers	Apache-2.0
Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause; BSD-3-Clause; MIT; CC-BY-2.5
Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause; MIT; public-domain; CC-BY-2.5
Guava GWT compatible libs		Apache-2.0; public-domain
Guava: Google Core Libraries for Java		public-domain; Apache-2.0
gwt-servlet		Apache-2.0; public-domain
gwt-user		Apache-2.0; public-domain
Hamcrest All		BSD-3-Clause-Clear
HAPI - 01 - Base Module		Multi-license: GPL-3.0-only OR MPL-1.1; Zlib
HAPI - Java HL7 API - Generated Structures v2.7		Multi-license: GPL-3.0-only OR MPL-1.1
HAPI - The Open Source Java HL7 Parser and API	Axel Biernat, Rodolfo Chavez, Ryan W. Gross, Nicolas Lefebvre, Jens Kristian Villadsen, Bryan Tripp, James Agnew, Christian Ohr, Niranjan Sharma, Alexei Guevara, Charles D. Fisher, Abraham Tio, Leslie Mann, Neal Acharya, Nico Vannieuwenhuyze	Multi-license: GPL-3.0-only OR MPL-1.1 OR MPL-2.0 OR Dual-license; Multi-license: GPL-3.0-only OR MPL-1.1
hazelcast		Apache-2.0; MIT
hazelcast-client		Apache-2.0
hazelcast-spring		Apache-2.0

HdrHistogram		Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause; public-domain
Hibernate Commons Annotations		LGPL-2.1-only; LGPL-3.0-only
Hibernate ORM - hibernate-core		LGPL-2.1-only; LGPL-2.1-or-later; LGPL-3.0-only
Hibernate ORM - hibernate-core		Multi-license: LGPL-2.1-only OR LGPL-2.1-or-later; LGPL-2.1-or-later; LGPL-3.0-only; EPL-2.0
Hibernate Validator Engine		Apache-2.0
Hibernate Validator Engine		Apache-2.0
HttpClient		Apache-2.0
HttpClient		Apache-2.0
HttpCore		Apache-2.0
ICU4J	Mark Davis, John Emmons, Doug Felt, Deborah Goldsmith, Steven Loomis, Markus Scherer, Peter Edberg, Yoshito Umaoka	Multi-license: BSD-3-Clause OR Unicode/ICU License; BSD-3-Clause; public-domain; SPL-1.0
itext	Bruno Lowagie, Paulo Soares, Mark Hall, Xavier Le Vourch	Multi-license: LGPL-3.0-only OR MPL-1.1 OR MPL-2.0; Multi-license: LGPL-3.0-only OR MPL-1.1; Multi-license: LGPL-2.0-or-later OR MPL-1.1; Multi-license: BSD-3-Clause OR LGPL-2.0-or-later OR MPL-1.1; Multi-license: LGPL-2.0-or-later OR MIT OR MPL-1.1; Multi-license: Apache-2.0 OR LGPL-2.0-or-later OR MPL-1.1; MPL-1.1; MIT; Multi-license: LGPL-3.0-only OR MIT OR MPL-2.0; Apache-2.0; LGPL-2.1-or-later; Multi-license: Apache-2.0 OR GPL-2.0-only OR LGPL-2.1-only; BSD-3-Clause; Multi-license: BSD-3-Clause OR SGI-B-2.0; SGI-B-2.0; CC-BY-2.5; LGPL-2.0-only; Adobe-2006; public-domain
Jackson datatype: jdk8		Apache-2.0
Jackson datatype: JSR310		Apache-2.0
Jackson-annotations		Apache-2.0
Jackson-core		Apache-2.0
jackson-databind		Apache-2.0

Jackson-dataformat-CSV		Apache-2.0
Jackson-dataformat-YAML		Apache-2.0
Jackson-dataformat-YAML		Apache-2.0
Jackson-datatype-Guava		Apache-2.0
JasperReports	Teodor Danciu, Lucian Chirita	LGPL-3.0-only; LGPL-3.0-or-later; GPL-3.0-only
JasperReports Font Extension		LGPL-3.0-only
JasperReports Functions		LGPL-3.0-only; LGPL-3.0-or-later
Java Annotation Indexer		Apache-2.0
Java Architecture for XML Binding		
Java Native Access	Timothy Wall, Matthias Bränding	Multi-license: Apache-2.0 OR LGPL-2.1-only
Java Persistence API, Version 2.1	Emmanuel Bernard, Hardy Ferentschik, Steve Ebersole	PHP-3.01; EPL-1.0
Java Servlet API	Rajiv Mordani, Shing Wai Chan	Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: Apache-2.0 OR CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; CDDL-1.1
JavaBeans(TM) Activation Framework		CDDL-1.1; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception
JavaMail API		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; CDDL-1.1
Javassist	Shigeru Chiba, Andrew Dinn, Kabir Khan, Scott Marlow	Multi-license: Apache-2.0 OR LGPL-2.1-only OR MPL-1.1 OR MPL-2.0; Multi-license: Apache-2.0 OR LGPL-2.1-or-later OR MPL-1.1; Multi-license: Apache-2.0 OR LGPL-2.1-only OR MPL-1.1

Javassist	Shigeru Chiba, Andrew Dinn, Kabir Khan, Scott Marlow	Multi-license: Apache-2.0 OR LGPL-2.1-only OR MPL-1.1 OR MPL-2.0; Multi-license: Apache-2.0 OR LGPL-2.1-only OR MPL-1.1; Multi-license: Apache-2.0 OR LGPL-2.1-or-later OR MPL-1.1
javax.annotation-api		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; CDDL-1.1
JAXB RI		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception OR public-domain
jaxb-api	Martin Grebac, Iaroslav Savvitskyi	Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; CDDL-1.1
jaxen	Ryan Gustafson, David Peterson, Mark Wilson, Jacob Kjome, Michael Brennan, Jason Hunter, Brett McLaughlin, Bradley S. Huffman, K. Ari Krupnikov, Paul R. Brown, Guoliang Cao, Jérôme Nègre, Eddie McGreal, Steen Lehmann, Ben McCann, Don Corley, Bob McWhirter, James Strachan, David Megginson, Erwin Bolwidt	http://jaxen.codehaus.org/licenses.html ; BSD-3-Clause; W3C
jCIFS		LGPL-2.1-only; LGPL-2.1-or-later; BSD-2-Clause; LGPL-2.0-or-later
jCSV		Apache-2.0
Jena		BSD-3-Clause; Apache-2.0; MS-PL

JGraphT - Core		Multi-license: EPL-1.0 OR LGPL-2.1-only; LGPL-2.1-or-later; LGPL-2.1-only
Joda-Time	Adrian Acala, Afif Ahmed, AlexMe951, Guy Allard, Brad Arndt, Mikey Austin, Oren Benjamin, Fredrik Borgh, Dave Brosius, Dan Cavallaro, Luc Claes, Emiliano Claria, Dan Cojocar, Evgeniy Devyatykh, dspitfire, Christopher Elkins, emopers, Michael Ernst, Jeroen van Erp, Gwyn Evans	Apache-2.0; public-domain
JSON Small and Fast Parser	Uriel Chemouni, Eitan Raviv	Apache-2.0
JSON.simple		Apache-2.0
jsoup		MIT
JUnit	David Saff, Kevin Cooney, Stefan Birkner, Marc Philipp	EPL-1.0
LatencyUtils		public-domain
Liquibase		Multi-license: Apache-2.0 OR BSD-3-Clause OR GPL-2.0-only; MIT; Apache-2.0
liquibase-slf4j		MIT
log4j		Apache-2.0
Lucene Common Analyzers		Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause; BSD-3-Clause; BSD-2-Clause; Multi-license: Apache-2.0 OR BSD-2-Clause
Lucene Core		Apache-2.0; BSD-3-Clause; MIT
Lucene QueryParsers		Apache-2.0

mariadb-java-client	Vladislav Vaintroub, Massimo Siani, Georg Richter	Multi-license: BSD-3-Clause OR LGPL-2.1-only OR LGPL-2.1-or-later; Multi-license: BSD-3-Clause OR LGPL-2.1-or-later
micrometer-core		Apache-2.0; Multi-license: Apache-2.0 OR MIT
micrometer-registry-new-relic		Apache-2.0
micrometer-registry-prometheus		Apache-2.0
Microsoft JDBC Driver for SQL Server		MIT; Apache-2.0; public-domain
mockito-core	Szczepan Faber, Brice Dutheil, Rafael Winterhalter, Tim van der Lippe	MIT; Multi-license: Apache-2.0 OR MIT; Apache-2.0; BSD-3-Clause
mockito-core	Szczepan Faber, Brice Dutheil, Rafael Winterhalter, Tim van der Lippe	MIT; Multi-license: Apache-2.0 OR MIT; Apache-2.0
Netty/All-in-One		Apache-2.0; BSD-3-Clause; MIT
OAuth2 for Spring Security		Apache-2.0
Objenesis		Apache-2.0
OGNL - Object Graph Navigation Library		Apache-2.0; BSD-3-Clause
Old JAXB Core		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception OR public-domain
org.apache.poi.xwpf.converter.pdf		MIT
OWASP Java HTML Sanitizer		Apache-2.0; BSD-3-Clause
Pirity User		Apache-2.0; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception
quartz		Apache-2.0
REngine Java interface to R		Multi-license: GPL-2.0-only OR LGPL-2.1-only
Rserve client back-end for REngine		Multi-license: GPL-2.0-only OR LGPL-2.1-only

Spring Beans		Apache-2.0
Spring Context		Apache-2.0
Spring Core		Apache-2.0; BSD-3-Clause
Spring HATEOAS	Oliver Gierke, Greg Turnquist	Apache-2.0
Spring JMS		Apache-2.0
Spring LDAP Core	Eric Dalquist, Marius Scurtescu, Tim Terry, Keith Barlow, Paul Harvey, Marvin S. Addison, Mattias Hellborg Arthursson, Ulrik Sandberg	Apache-2.0
Spring Messaging		Apache-2.0
Spring Object/Relational Mapping		Apache-2.0
Spring TestContext Framework		Apache-2.0
Spring Transaction		Apache-2.0
Spring Web		Apache-2.0; public-domain
Spring Web		Apache-2.0; public-domain
Spring Web MVC		Apache-2.0
spring-boot-starter-thymeleaf		Apache-2.0
spring-boot-starter-web		Apache-2.0
spring-security-config	Rob Winch, Joe Grandja	Apache-2.0
spring-security-core	Rob Winch, Joe Grandja	Apache-2.0
spring-security-ldap	Rob Winch, Joe Grandja	Apache-2.0
spring-security-oauth2-client	Rob Winch, Joe Grandja	Apache-2.0
spring-security-oauth2-jose	Rob Winch, Joe Grandja	Apache-2.0
spring-security-saml2-service-provider	Rob Winch, Joe Grandja	Apache-2.0
spring-security-web	Rob Winch, Joe Grandja	Apache-2.0

springfox-core	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-swagger-ui	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-swagger2	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
swagger-annotations		Apache-2.0
swagger-models		Apache-2.0
thymeleaf	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0
thymeleaf-spring5	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0
unbescape		Apache-2.0
vaadin-server		Apache-2.0
vaadin-spring		Apache-2.0
vaadin-themes		Apache-2.0
VT Password Libraries	Daniel Fisher, Marvin Addison	Multi-license: Apache-2.0 OR LGPL-2.1-only OR LGPL-3.0-only
Xalan Java		W3C; Apache-2.0
XBean :: Spring		Apache-2.0
Xerces2-j		Apache-2.0; W3C
ZXing Core		Apache-2.0
Deep Dependencies		
Title	Authors	Declared License
ActiveMQ :: Broker		Apache-2.0
ActiveMQ :: Client		Apache-2.0
ActiveMQ :: Generic JMS Pool		Apache-2.0
ActiveMQ :: Openwire Legacy Support		Apache-2.0
ActiveMQ :: Pool		Apache-2.0
ActiveMQ :: Spring		Apache-2.0
ActiveMQ Protocol Buffers Implementation and Compiler		Apache-2.0
antlr		Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause

AOP alliance		public-domain
Apache Ant + ANTLR		Apache-2.0
Apache Ant + JUnit		Apache-2.0
Apache Ant Core		Apache-2.0
Apache Ant Launcher		Apache-2.0
Apache Commons Codec	Christopher O'Brien, Martin Redington, Jeffery Dever, Steve Zimmermann, Benjamin Walstrum, Oleg Kalnichevski, Dave Dribin, Alex Karasulu, Matthew Inger, Jochen Wiedmann, Sebastian Bazley, Matthew Pocock, Colm Rice, Henri Yandell, Tim O'Brien, Scott Sanders, Rodney Waldhoff, Daniel Rall, Jon S. Stevens, Gary Gregory	Apache-2.0
Apache Commons Collections	Rafael U. C. Afonso, Max Rydahl Andersen, Avalon, Federico Barbieri, Jeffrey Barnes, Nicola Ken Barozzi, Arron Bates, Sebastian Bazley, Benjamin Bentmann, Ola Berg, Sam Berlin, Christopher Berry, Nathan Beyer, Rune Peter Bjørnstad, Janek Bogucki, Maarten Brak, Dave Bryson, Chuck Burdick, Julien Buret, Josh Cain	Apache-2.0

Apache Commons Collections	Rafael U. C. Afonso, Max Rydahl Andersen, Avalon, Federico Barbieri, Jeffrey Barnes, Nicola Ken Barozzi, Arron Bates, Sebastian Bazley, Benjamin Bentmann, Ola Berg, Sam Berlin, Christopher Berry, Nathan Beyer, Rune Peter Bjørnstad, Janek Bogucki, Maarten Brak, Dave Bryson, Chuck Burdick, Julien Buret, Josh Cain	Apache-2.0
Apache Commons Compress	Wolfgang Glas, Christian Kohlschütter, Bear Giles, Michael Kuss, Lasse Collin, John Kodis, BELUGA BEHR, Simon Spero, Michael Hausegger, Torsten Curdt, Stefan Bodewig, Sebastian Bazley, Christian Grobmeier, Julius Davies, Damjan Jovanovic, Emmanuel Bourg, Gary Gregory, Rob Tompkins, Peter Alfred Lee	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause
Apache Commons Compress	Wolfgang Glas, Christian Kohlschütter, Bear Giles, Michael Kuss, Lasse Collin, John Kodis, BELUGA BEHR, Simon Spero, Michael Hausegger, Torsten Curdt, Stefan Bodewig, Sebastian Bazley, Christian Grobmeier, Julius Davies, Damjan Jovanovic, Emmanuel Bourg, Gary Gregory, Rob Tompkins, Peter Alfred Lee	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause
Apache Commons IO	Rahul Akolkar, Jason Anderson, Nathan Beyer, Emmanuel Bourg, Chris Eldredge, Magnus Grimsell, Jim Harrington, Thomas Ledoux, Andy Lehane, Marcelo Liberato, Alban Peignier, Adam Retter, Ian Springer, Dominik Stadler, Masato Tezuka, James Urie, Frank W. Zammetti, Scott Sanders, dlon Gillard, Nicola Ken Barozzi	Apache-2.0
Apache Commons Lang	C. Scott Ananian, Chris Audley, Stephane Bailliez, Michael Becke, Benjamin Bentmann, Ola Berg, Nathan Beyer, Stefan Bodewig, Janek Bogucki, Mike Bowler, Sean Brown, Alexander Day Chaffee, Al Chou, Greg Coladonato, Maarten Coene, Justin Couch, Michael Davey, Norm Deane, Morgan Delagrance, Ringo De Smet	Apache-2.0
Apache Commons Lang	C. Scott Ananian, Chris Audley, Stephane Bailliez, Michael Becke, Benjamin Bentmann, Ola Berg, Nathan Beyer, Stefan Bodewig, Janek Bogucki, Mike	Apache-2.0

	Bowler, Sean Brown, Alexander Day Chaffee, Al Chou, Greg Coladonato, Maarten Coene, Justin Couch, Michael Davey, Norm Deane, Morgan Delagrange, Ringo De Smet	
Apache Commons Math	Eldar Agalarov, Tim Allison, C. Scott Ananian, Mark Anderson, Peter Andrews, RÅ@mi Arntzen, Matt Adereth, Jared Becksfort, Michael Bjorkegren, Brian Bloniarz, John Bollinger, Cyril Briquet, Dave Brosius, Dan Checkoway, Anders Conbere, Charles Cooper, Paul Cowan, Benjamin Croizet, Larry Diamond, Aleksei Dievskii	Multi-license: Apache-2.0 OR BSD-3-Clause; BSD-2-Clause; BSD-3-Clause; Multi-license: Apache-2.0 OR BSD-2-Clause; Apache-2.0
Apache Commons Net	Henrik Sorensen, Jeff Nadler, William Noto, Stephane ESTE-GRACIAS, Dan Armbrust, Yuval Kashtan, Joseph Hindsley, Rob Hasselbaum, Mario Ivankovits, Naz Irizarry, Tapan Karecha, Jason Mathews, Winston Ojeda, Ted Wise, Bogdan Drozdowski, Jeffrey D. Brekke, Steve Cohen, Bruno D'Avanzo, Daniel F. Savarese, Rory Winston	Apache-2.0; Multi-license: Apache-2.0 OR NTP
Apache Commons Pool	Morgan Delagrange, Geir Magnusson, Craig McClanahan, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Robert Burrell Donkin, Sandy McArthur, Simone Tripodi, Gary Gregory, Matt Sicker	Apache-2.0
Apache Commons Pool	Morgan Delagrange, Geir Magnusson, Craig McClanahan, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Robert Burrell Donkin, Sandy McArthur, Simone Tripodi, Gary Gregory, Matt Sicker	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0; public-domain
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian	Apache-2.0

	McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0; Multi-license: Apache-2.0 OR CC-BY-2.5
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll,	Apache-2.0; MIT

	Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0; MIT
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian	Apache-2.0

	McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0

Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0

Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udoenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	Apache-2.0

	Ganza, Arjun Nayyar, Mark Chu-Carroll, Mark Turansky, Jean-Louis Berliet, Graham Miller, Marc Palmer	
Apache JempBox		Apache-2.0
Apache POI		Apache-2.0
Apache Preflight		Apache-2.0
Apache Tika core		Apache-2.0
Apache XML Graphics Commons		Apache-2.0
Apache XML Security for Java	Sean Mullan, Berin Lautenbach, Axl Mattheus, Erwin van der Koog, Raul Benito Garcia, Vishal Mahajan, Milan Tomic, Colm O hEigeartaigh	Apache-2.0; W3C
Apache XmpBox		Apache-2.0
asm	Eric Bruneton, Eugene Kuleshov, Remi Forax	Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause
attoparser		Apache-2.0
Avalon Framework Implementation		Apache-2.0
batik-anim		Apache-2.0
batik-constants		Apache-2.0
batik-css		Apache-2.0
batik-ext		Apache-2.0; W3C
batik-gvt		Apache-2.0
batik-i18n		Apache-2.0
batik-parser		Apache-2.0
batik-script		Apache-2.0
batik-shared-resources		Apache-2.0
batik-svg-dom		Apache-2.0; W3C
BeanUtils	Robert Burrell Donkin, dlon Gillard, Craig McClanahan, Geir Magnusson Jr., Scott Sanders, James Strachan, Rodney Waldhoff, Martin van den Bemt	Apache-2.0
Bouncy Castle ASN.1 Extension and Utility APIs		Bouncy Castle Licence
Bouncy Castle PKIX, CMS, EAC, TSP, PKCS, OCSP, CMP, and CRMF APIs		Bouncy Castle Licence; public-domain
Byte Buddy (without dependencies)		Apache-2.0; BSD-3-Clause
c3p0		Multi-license: LGPL-2.1-only OR EPL-1.0; LGPL-2.1-only
Castor CORE - Core code/functionality		Apache-2.0; BSD-3-Clause; unknown
Castor XML - core		Apache-2.0; BSD-3-Clause; unknown
CGLIB		Apache-2.0; BSD-3-Clause
Checker Qual	Michael Ernst, Suzanne Millstein	MIT; SPL-1.0

ClassMate	Tatu Saloranta, Brian Langel	Apache-2.0; unknown
ClassMate	Tatu Saloranta, Brian Langel	Apache-2.0; unknown
Collections	Max Rydahl Andersen, Federico Barbieri, Arron Bates, Nicola Ken Barozzi, Ola Berg, Christopher Berry, Janek Bogucki, Chuck Burdick, Dave Bryson, Julien Buret, Jonathan Carlson, Ram Chidambaram, Peter Donald, Steve Downey, Rich Dougherty, Stefano Fornari, Andrew Freeman, Gerhard Froehlich, Paul Jack, Eric Johnson	Apache-2.0
Commons Digester		Apache-2.0
Commons Digester		Apache-2.0
Commons Pool	Morgan Delagrance, Geir Magnusson, Craig McClanahan, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Robert Burrell Donkin, Sandy McArthur, Phil Steitz	Apache-2.0
Commons Pool	Morgan Delagrance, Geir Magnusson, Craig McClanahan, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Robert Burrell Donkin, Sandy McArthur, Phil Steitz	Apache-2.0

Cryptacular Library	Daniel Fisher, Marvin S. Addison	Multi-license: Apache-2.0 OR LGPL-3.0-only
curvesapi	stormdollar, Dustin Spicuzza	Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause; Apache-2.0
dec		MIT
dom4j		Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause
Eclipse Compiler for Java(TM)		Multi-license: EPL-1.0 OR EPL-2.0; EPL-2.0; EPL-1.0
Elemental2 Promise		Apache-2.0
error-prone annotations		Apache-2.0
Extended StAX API	Roman Grigoriadi, Zheng Jun Li	Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception
fastinfoset		Apache-2.0; Multi-license: Apache-2.0 OR BSD-2-Clause; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception
Flute		W3C; Multi-license: GPL-2.0-only OR W3C
fr.opensagres.poi.xwpf.converter.core		MIT
fr.opensagres.poi.xwpf.converter.pdf		MIT
fr.opensagres.poi.xwpf.converter.xhtml		MIT
fr.opensagres.xdocreport.converter		MIT
fr.opensagres.xdocreport.core		MIT; Multi-license: Apache-2.0 OR MIT
fr.opensagres.xdocreport.itext.extension		MIT
GenTyRef		Apache-2.0
Google Guice - Core Library		Apache-2.0
Google Guice - Extensions - AssistedInject		Apache-2.0
Guava InternalFutureFailureAccess and InternalFutures		Apache-2.0
Guava ListenableFuture only		Apache-2.0

Hamcrest Core		
hawtbuf		Apache-2.0
Hibernate Commons Annotations		Multi-license: LGPL-2.1-only OR LGPL-2.1-or-later; LGPL-2.1-or-later
HikariCP		Apache-2.0
HttpClient		Apache-2.0
HttpCore		Apache-2.0
ICU4J	Mark Davis, John Emmons, Doug Felt, Deborah Goldsmith, Steven Loomis, Markus Scherer, Peter Edberg, Yoshito Umaoka	ICU; MIT; GPL-3.0-only
IRI		BSD-3-Clause
istack common utility code runtime		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception
J2EE Management 1.1		Apache-2.0
J2ObjC Annotations		Apache-2.0
Jackson BOM		Apache-2.0
Jackson-annotations		Apache-2.0
Jackson-core		Apache-2.0
jackson-databind		Apache-2.0
jakarta-regexp		Apache-2.0
Java Advanced Imaging Image I/O Tools API core (standalone)	John Hewson, Peter Hull, Mark Carroll, Robin Stevens, Yannick De Turck, Butch Howard	Multi-license: BSD-3-Clause OR GPL-2.0-only; Custom License File
Java Annotation Indexer		Apache-2.0
Java Servlet API	Rajiv Mordani, Shing Wai Chan	Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: Apache-2.0 OR CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; CDDL-1.1

Java Transaction API		Multi-license: GPL-2.0-with-classpath-exception OR Common Development and Distribution License; GPL-2.0-with-classpath-exception
java-support		Multi-license: Apache-2.0 OR BSD-3-Clause; Apache-2.0
JavaBeans Activation Framework API jar		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; CDDL-1.1
JavaBeans(TM) Activation Framework		CDDL-1.1
JavaMail API		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; GPL-2.0-only; Multi-license: GPL-2.0-only OR GPL-2.0-with-classpath-exception; CDDL-1.1
JavaServlet(TM) Specification		Apache-2.0
javax.inject		Apache-2.0
javax.persistence-api		EPL-1.0
JAXB RI		Multi-license: CDDL-1.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception OR GPL-3.0-only; Multi-license: CDDL-1.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception OR Public-Domain
JAXB Runtime		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception OR public-domain
JBoss Logging 3		Apache-2.0
JCIP Annotations under Apache License		Apache-2.0
JCL 1.2 implemented over SLF4J		Apache-2.0
JCL 1.2 implemented over SLF4J		MIT; Apache-2.0

JCommander		Apache-2.0
JCommon		LGPL-3.0-only; LGPL-2.1-or-later; Multi-license: GPL-2.0-only OR LGPL-2.1-or-later; GPL-2.0-only; LGPL-2.1-only
JDOM	Jason Hunter, Rolf Lear	Apache-2.0; BSD-3-Clause
JFreeChart		LGPL-3.0-only; LGPL-2.1-or-later; BSD-3-Clause
JLine	Scott Gray, Hiram Chirino, Jason Dillon, Guillaume Nodet, Marc Prud'hommeaux	Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause
JMS 1.1		Apache-2.0
Joda-Time	Adrian Acala, Afif Ahmed, AlexMe951, Guy Allard, Brad Arndt, Mikey Austin, Oren Benjamin, Fredrik Borgh, Dave Brosius, Dan Cavallaro, Luc Claes, Emiliano Claria, Dan Cojocar, Evgeniy Devyatkykh, dspitfire, Christopher Elkins, emopers, Michael Ernst, Jeroen van Erp, Gwyn Evans	Apache-2.0; public-domain
JsInterop Base		Apache-2.0
jsinterop-annotations		Apache-2.0
jsinterop-annotations		Apache-2.0
jsoup		MIT
JTA 1.1		Apache-2.0
JTidy		Java HTML Tidy License; public-domain; W3C
JUL to SLF4J bridge		MIT
JUL to SLF4J bridge		MIT
JUnit Jupiter API	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0; unknown
JUnit Jupiter Engine	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0; unknown
JUnit Platform Commons	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0; unknown
JUnit Platform Engine API	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0; unknown
JUnit Platform Launcher	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0; unknown
Log4j Implemented Over SLF4J		Apache-2.0
Lucene Queries		Apache-2.0

Lucene Sandbox		Apache-2.0
MapStruct Core	Gunnar Morling, Filip Hrisafov	Apache-2.0
mchange-commons-java		Multi-license: EPL-1.0 OR LGPL-2.1-only; LGPL-2.1-only; Multi-license: GPL-2.0-only OR LGPL-2.1-only
Metrics Core		public-domain
Nimbus Content Type		Apache-2.0
Nimbus JOSE+JWT		Apache-2.0
Nimbus LangTag		Apache-2.0
OAuth 2.0 SDK with OpenID Connect extensions		Apache-2.0
Objenesis		Apache-2.0
OGNL - Object Graph Navigation Library		Apache-2.0; BSD-3-Clause
Old JAXB Core		Multi-license: CDDL-1.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception OR GPL-3.0-only; Multi-license: CDDL-1.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception OR Public-Domain
OOXML schemas		Apache-2.0
OpenSAML :: Core		Apache-2.0
OpenSAML :: Messaging API		Apache-2.0
OpenSAML :: Profile API		Apache-2.0
OpenSAML :: SAML Provider API		Apache-2.0
OpenSAML :: SAML Provider Implementations		Apache-2.0
OpenSAML :: Security API		Apache-2.0
OpenSAML :: Security Implementation		Apache-2.0
OpenSAML :: SOAP Provider API		Apache-2.0; W3C
OpenSAML :: SOAP Provider Implementations		Apache-2.0
OpenSAML :: Storage API		Apache-2.0
OpenSAML :: XML Security API		W3C; Apache-2.0
OpenSAML :: XML Security Implementation		Apache-2.0

org.apache.poi.xwpf.converter.core		MIT
org.opentest4j:opentest4j	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen	Apache-2.0
parso	Petr Tsurinov, Igor Printsev	Apache-2.0
PDFBox JBIG2 ImageIO plugin	Andreas Lehmkuhler, Adam Nichols, Ben Litchfield, Brian Carrier, Daniel Wilson, Guillaume Bailleul, Jürg Henne, Sebastian Holder, John Hewson, Jeremias Maerki, Kevin Jackson, Johannes Koch, Carolin Köhler, Matthias Mayer, Maruan Sayhoun, Phillipp Koch, Tim Allison, Thomas Chojecki, Timo Boehme, Tilman Hausherr	Apache-2.0
picocli - a mighty tiny Command Line Interface		Apache-2.0
Prometheus Java Simpleclient		Apache-2.0; public-domain
Prometheus Java Simpleclient Common		Apache-2.0
QDox	Mike Royle, Peter Donald, James Strachan, Nick Pomfret, Chris Stevenson, Ben Hogan, Laurent Etiemble, Shawn Chain, Brian Slesinsky, James Lee, Eric Redmond, Joe Walnes, Aslak Hellesoy, Paul Hamman, Mike Williams, Mauro Talevi, Robert Scholte	Apache-2.0
SentimentAnalysisParser		Apache-2.0
Simple API for CSS		W3C
SLF4J API Module		MIT
SLF4J API Module		MIT
SLF4J API Module		MIT
SLF4J API Module		MIT
SnakeYAML	Andrey Somov, Alexander Maslov, Jordan Angold	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR EPL-1.0 OR GPL-2.0-or-later OR LGPL-2.1-or-later

SnakeYAML	Andrey Somov, Alexander Maslov, Jordan Angold	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR EPL-1.0 OR GPL-2.0-or-later OR LGPL-2.1-or-later
SparseBitSet		Apache-2.0
Spring AOP		Apache-2.0
Spring AOP		Apache-2.0
Spring AOP		Apache-2.0
Spring Beans		Apache-2.0
Spring Beans		Apache-2.0
Spring Boot Tomcat Starter		Apache-2.0
Spring Commons Logging Bridge		Apache-2.0
Spring Commons Logging Bridge		Apache-2.0
Spring Context		Apache-2.0
Spring Context		Apache-2.0
Spring Core		Apache-2.0; BSD-3-Clause
Spring Core		Apache-2.0; BSD-3-Clause
Spring Expression Language (SpEL)		Apache-2.0
Spring Expression Language (SpEL)		Apache-2.0
Spring Expression Language (SpEL)		Apache-2.0
Spring JDBC		Apache-2.0
Spring Plugin - Metadata Extension		Apache-2.0
Spring Plugin Core		Apache-2.0
Spring Transaction		Apache-2.0
Spring Transaction		Apache-2.0
Spring Web		Apache-2.0
Spring Web MVC		Apache-2.0
Spring Web MVC		Apache-2.0
spring-boot		Apache-2.0
spring-boot-autoconfigure		Apache-2.0
spring-boot-starter		Apache-2.0
spring-boot-starter-logging		Apache-2.0
spring-ldap-core	Eric Dalquist, Marius Scurtescu, Tim Terry, Keith Barlow, Paul Harvey, Marvin S. Addison, Rob Winch, Mattias Hellborg Arthursson, Ulrik Sandberg	Apache-2.0

spring-security-crypto	Rob Winch, Joe Grandja	Apache-2.0; MIT
spring-security-oauth2-core	Rob Winch, Joe Grandja	Apache-2.0
springfox-schema	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-spi	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-spring-web	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-swagger-common	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
StAX API	Aleksander Slominski, Chris Fry	Apache-2.0; public-domain
Stax2 API		Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause
telemetry		Apache-2.0
telemetry-core		Apache-2.0
testng	Cedric Beust, Julien Herr, Krishnan Mahadevan	Apache-2.0
The Netty Project		Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause; Multi-license: Apache-2.0 OR MIT; BSD-3-Clause; Multi-license: LGPL-2.1-only OR MIT; LGPL-2.1-only; MIT; Public-Domain
thymeleaf	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0; W3C; IOS
thymeleaf-layout-dialect	Justin Munn, Oliver Niekrenz	Apache-2.0
thymeleaf-spring4	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0; W3C
tomcat-annotations-api		Apache-2.0

tomcat-embed-core		Apache-2.0; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; CDDL-1.1
tomcat-embed-el		Apache-2.0
tomcat-embed-websocket		Apache-2.0
TXW2 Runtime		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception OR public-domain
unbescape		Apache-2.0
Vaadin Sass Compiler		Apache-2.0; Multi-license: Apache-2.0 OR W3C
vaadin-shared		Apache-2.0
VT Crypt Library	Daniel Fisher, Marvin Addison	Multi-license: Apache-2.0 OR LGPL-2.1-only OR LGPL-3.0-only
VT Dictionary Libraries	Daniel Fisher, Marvin Addison	Multi-license: Apache-2.0 OR LGPL-2.1-only OR LGPL-3.0-only
Woodstox		Apache-2.0
Woodstox		Apache-2.0; LGPL-2.1-only
Xalan Java Serializer		Apache-2.0; W3C
Xerces2-j		Apache-1.1; W3C; BSD-2-Clause
XHTMLRenderer core-renderer		LGPL-2.1-only; Public-Domain
XML Commons External Components XML APIs Extensions		Apache-2.0
xml-apis		W3C; Apache-2.0; Multi-license: Apache-2.0 OR The SAX License; public-domain
XmiBeans	Cezar Andrei, Radu Preotiu, Wing Yew Poon, Jacob Danner, POI Team	Apache-2.0; Apache-1.1

Learn more:

www.agilent.com/chem/agilent Slims

U.S. and Canada

1-800-227-9770 agilent_inquiries@agilent.com

Europe

info_agilent@agilent.com

Asia Pacific

inquiry_lsca@agilent.com

This information is subject to change without notice.

DE44456.1092476852

© Agilent Technologies, Inc. 2021

Published in Europe, December 7, 2022

5994-4171EN

