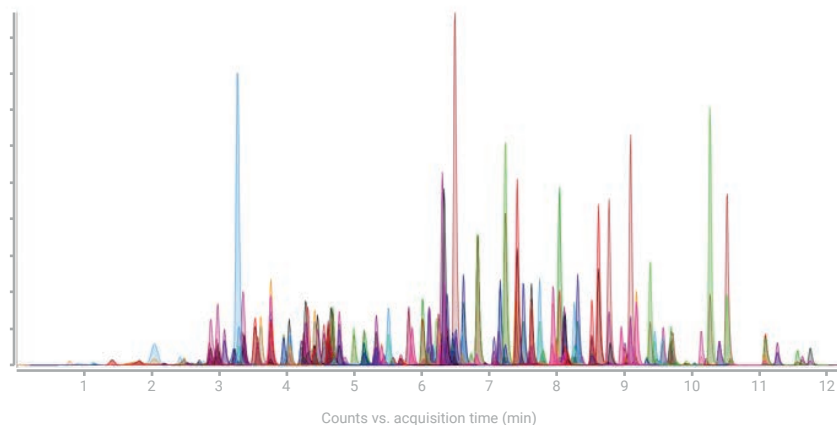


Multi-class, Multi-residue Veterinary Drug Analysis in Food Matrices

Everything you need to analyze 210 veterinary drugs with confidence

Agilent multi-class veterinary drug products help you simplify laboratory operations and combine multiple food analyses into one easy-to-follow protocol. So you can get up and running fast, maintain a well-trained staff, and maximize your productivity. Our portfolio includes:

- Optimized method, validated per AOAC, EU, and U.S. FDA guidelines
- Demonstrated method performance, reproducibility, and transferability
- Consistent sample preparation workflow, dMRM methods, and database
- Predefined consumables and 25-compound system suitability standard
- Comprehensive workflow guide complemented by onsite and online training
- Compatibility with Agilent 6470 and 6495C triple quadrupole LC/MS systems



Meet challenging AOAC, EU, and U.S. FDA screening requirements. This chromatogram represents the analysis of 210 veterinary drugs in chicken muscle.

Easy ordering information

Want to add the items under each category to your My Favorites¹ list at the Agilent online store? Simply click the header links below. Then, enter the quantities for the products you need. Your list of items will remain under My Favorites for you to use with future orders.

Description	Part Number
HPLC columns and supplies	
InfinityLab Poroshell 120 EC-C18, 3.0 x 100 mm, 2.7 µm, 1000 bar	695575-302
InfinityLab Poroshell 120 EC-C18, 3.0 x 5 mm, 2.7 µm, guard	823750-911
1290 Infinity II inline filter (0.3 µm)	5067-6189
Quick connect assembly, 0.12 x 105 mm, for connection at column inlet on LC system	5067-5957
Quick connect assembly, 0.17 x 105 mm, for connection at column inlet on LC system	5067-6166
Quick Turn fitting, for connection between column outlet and MS detector	5067-5966
Quick Turn capillary, 0.12 x 280 mm, for column outlet	5500-1191
Sample preparation	
Captiva EMR-Lipid, 3 mL cartridges, with 300 mg sorbent mass, 50/pk	5190-1003
Captiva EMR-Lipid, 6 mL cartridges	5190-1004 ²
Ceramic homogenizers, 50 mL tubes, 100/pk	5982-9313
Ceramic homogenizers, 15 mL tubes, 100/pk (optional)	5982-9312
50 mL plastic centrifuge tubes	5610-2049
Agilent positive pressure manifold 48 processor	5191-4101 ³
3 mL SPE cartridge rack, PPM-48	5191-4103 ³
6 mL SPE cartridge rack, PPM-48	5191-4104 ^{2,3}
Collection rack, 13 x 100 mg tubes, PPM-48	5191-4107 ³
Collection rack, 16 x 100 mm tubes, PPM-48	5191-4108 ³
Vials and caps	
Vial, amber, write-on spot, certified, 2 mL, 100/pk; vial size: 12 x 32 mm (12 mm cap)	5182-0716
Vial, high recovery, amber, 1.5 mL with 30 µL reservoir, 100/pk; vial size: 12 x 32 mm (12 mm cap)	5183-2073
Screw cap, bonded, blue, PFTE/red silicone septa, 100/pk; cap size: 12 mm	5190-7024
Screw vial, fixed insert, amber 12 x 32 mm	5188-6592
Vial rack, 12 mm vials, 5/pk	9301-0722

Description	Part Number
Solvents and reagents	
InfinityLab ultrapure LC/MS acetonitrile	5191-4496
InfinityLab ultrapure LC/MS methanol	5191-4497
InfinityLab ultrapure LC/MS water	5191-4498
5 M ammonium formate solution	G1946-85021
MS solution, formic acid, 99.5% purity, 10 mL	US-700002341
Formic acid reagent grade, 99.5% purity, 5 mL	G2453-85060
Chemical standards	
Vet drug system suitability mix: 25 component	5799-0015

Use the Agilent Veterinary Drug Suitability Mix as a system and method performance check for your instrument, column, and consumables. These mixes can be run daily, weekly, or monthly to maintain confidence in your results and identify issues early.

For more individual vet drug standards, visit Agilent [chemical standards](#).

¹ If this is the first time you are using "My Favorites," you will be asked to enter your email address for account verification. If you have an existing Agilent account, you will be able to log in. If you don't have a registered Agilent account, you will need to [register](#) for one. This feature is valid only in regions that are e-commerce enabled. All items can also be ordered through your regular sales and distributor channels.

² Either 3 or 6 mL can be used. Match cartridge size to cartridge rack size.

³ One-time purchase.

Effectively detect a range of veterinary drugs, metabolites, and transformation products in foods of animal origin.

Visit www.agilent.com/chem/vetdrugslctq

DE.4448958333

This information is subject to change without notice.