

Agilent BioTek Dual-Reagent Injector Module Chemical Compatibility



The Agilent BioTek dual-reagent injector module is available for use with the Agilent BioTek Synergy Neo2 hybrid multimode reader, Synergy H1 and HTX multimode readers, Cytation 1/5/7/9 cell imaging multimode readers, and Cytation C10 confocal imaging reader.

This technical overview describes the materials used in the module and information to assist in determining their compatibility with common chemicals.

Table 1. Locations of materials used in the Agilent BioTek dual-reagent injector module.

	Material	Location
1.	316 stainless steel	Inlet tube, injector tip
2.	Polyphenylene sulfide (PPS)	Valve body
3.	Ethylene propylene (EPDM)	Valve diaphragm
4.	Polytetrafluoroethylene (PTFE)	Syringe piston, all tubing
5.	Borosilicate glass	Syringe cylinder
6.	Chlorinated polyvinyl chloride (CPVC)	Injector block

Table 2. Chemical compatibility ratings.

Chemical	Material					
	1	2	3	4	5	6
Acetonitrile	A	A	D	A	A	D
Chloroform	A	A	D	A	A	D
Detergents (1%)	A	A	A	A	A	A
DMSO	A	A	B	A	A	D
Ethanol (70%)	A	A	A	A	A	B
Ethylene oxide	B	D	C	A	A	C
Formaldehyde (40%)	A	A	A	A	A	A
Hydrochloric acid (20%)	D	D	A	A	A	A
Hydrogen peroxide (10%)	B	A	A	A	A	A
Isopropyl alcohol (70%)	A	A	A	A	A	B
Methanol (70%)	A	A	A	A	A	A
Sodium hydroxide (20%)	B	A	B	A	A	A
Sodium hypochlorite (< 20%)	C	A	B	A	A	A
Sodium hypochlorite (0.5%)	B	A	B	A	A	A
Sulfuric acid (< 10%)	B	A	A	A	A	A
Virkon (10%)	A	ND	A	ND	A	A

Key: A = no effect; B = slight effect; C = moderate effect; D = severe effect; ND = no data

Note: Table 2 does not include all chemicals that can cause severe damage to one or more materials used in the system. Continuous contact with harsh chemicals is not recommended. It is best to rinse the system with de-ionized water after contact with any strong acid, base, or solvent. Table 2 does not list all possible chemicals and concentrations and should be used only as a guide. Some chemicals are known to cause severe damage to materials used in the injection system, including high concentrations of hydrochloric acid, undiluted sodium hypochlorite, and high concentrations of sulfuric acid. If there are questions about the compatibility of any chemicals not described, please contact Agilent.

www.agilent.com/lifesciences/biotek

DE-012150

This information is subject to change without notice.

© Agilent Technologies, Inc. 2026
Published in the USA, March, 2026
5994-6861EN

PN 1321028 Rev G