



CHROMATOGRAPHY

Chemical Compatibility with the NGC Chromatography System

The following table shows the chemicals that are compatible with the NGC Chromatography System.

- All materials used in the manufacture of the NGC System have met the required quality specifications and are safe for use with these chemicals
- The wettable parts of the NGC System are composed of polyetheretherketone (PEEK) plastic, high density polyethylene, and quartz glass. Materials used in the wetted parts of the NGC System have known chemical compatibilities, which are suitable for most common aqueous liquid chromatography applications and some organic liquid chromatography applications (such as reverse phase)
- The following chemicals can be used with the NGC System but long-term storage should be avoided:
 - Stronger acids
 - Aldehydes
 - Esters
 - Aliphatic hydrocarbons
 - 100% acetonitrile
- **Warning:** The following solvents will damage the NGC and are not recommended for even short-term use with the system:
 - Aromatic and halogenated hydrocarbons
 - Ketones
 - Strong oxidizing agents

Chemicals That Are Compatible with the NGC Chromatography System

Acids and Bases	Salts	Alcohols	Organics	Buffers	Detergents and Other
Acetic acid	6 M guanidine hydrochloride	100% ethanol	100% acetone	Bicine	0.1% bleach
Citric acid	Potassium chloride	40% ethylene glycol	100% acetonitrile	Bis-Tris	β-mercaptoethanol (β-ME)
Dimethyl arsenic acid (cacodylate)	Potassium dihydrogen orthophosphate	100% isopropanol	Diethanolamine	Bis-Tris hydrochloride	10% CHAPS
Formic acid	Potassium dihydrogen phosphate	100% methanol	Ethanolamine	Bis-Tris propane	Dithiothreitol (DTT)
2 M hydrochloric acid	Potassium hydrogen phosphate		100% methylene chloride	Glycine hydrochloride	50% EDTA
2 M sodium hydroxide	Potassium hydrogen phthalate		N-methyl piperazine	HEPES	40% glycerol
Succinic acid	Potassium phosphate		N-methyl piperazine dihydrochloride	MES monohydrate	30% hydrogen peroxide
1 M sulfuric acid	Saline sodium citrate		Organic amines	MOPS	2% lysozyme
1% trifluoroacetic acid	Sodium acetate trihydrate		Piperazine	PIPES	N-dodecyl-β-D-maltoside
	Sodium bicarbonate		Triethanolamine	Sodium formate	N-octyl-β-D-glucopyranoside
	10% sodium bromide		Triethanolamine hydrochloride	Sodium HEPES	10% SDS
	Sodium carbonate			Sodium MES	Sodium barbitone
	4 M sodium chloride			Sodium TAPS	8 M Tris (2-carboxyethyl) phosphine (TCEP)
	Sodium citrate			Sodium tetraborate	2% Triton X-100
	Sodium dihydrogen orthophosphate			TAPS	8 M urea
	Sodium dihydrogen phosphate			TES	
	Sodium hydrogen orthophosphate			Tricine	
	Sodium hydrogen phosphate			Tris base	
	Sodium hydrogen phthalate			Tris hydrochloride	
	Sodium phosphate dibasic heptahydrate			Trisodium citrate dihydrate	
	Sodium phosphate monobasic monohydrate				

Visit [bio-rad.com/NGCcompatibility](https://www.bio-rad.com/NGCcompatibility) for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.
All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site [bio-rad.com](https://www.bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00
Germany 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600
Japan 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30
Poland 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580
Sweden 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

