



Chemical Compatibility

	Material			Type
	RC membrane	Housing PP	Claristep® RC	
Filter Membrane	RC			RC
Housing Material		PP		PP
Sterilization				
Ethylene oxide	++	++	++	
Gamma irradiation	++	-	-	
Autoclaving 121°C, 30 min	++	++	++	
Solvents				
Acetone	++	++	++	
Acetonitrile	++	++	++	
Gasoline	++	++	++	
Benzene	++	++	++	
Benzyl alcohol	++	+	+	
n-Butyl acetate	++	++	++	
n-Butanol	++	++	++	
Cellosolve	++	++	++	
Chloroform	++	++	++	
Cyclohexane	++	+	+	
Cyclohexanone	++	+	+	
Diethylacetamide	++	++	++	
Diethyl ether	++	++	++	
Dimethyl formamide	+	++	+	
Dimethylsulfoxide	++	++	++	
Dioxane	++	++	++	
Ethanol, 98%	++	++	++	
Ethyl acetate	++	+	+	
Ethylene glycol	++	++	++	
Formamide	+	++	+	
Glycerin	++	++	++	
n-Heptane	++	+	+	
n-Hexane	++	+	+	
Isobutanol	++	++	++	
Isopropanol	++	++	++	
Isopropyl acetate	++	++	++	
Methanol, 98%	++	++	++	
Methyl acetate	++	+	+	
Methylene chloride	++	++	++	
Methyl ethyl ketone	++	+	+	
Methyl isobutyl ketone	++	+	+	
Monochlorobenzene	++	+	+	
Nitrobenzene	++	+	+	
n-Pentane	++	+	+	
Perchloroethylene	++	+	+	

	Material			Type
	RC membrane	Housing PP	Claristep® RC	
Filter Membrane	RC			RC
Housing Material		PP		PP
Solvents (continued)				
Pyridine	++	++	++	
Carbon tetrachloride	++	+	+	
Tetrahydrofuran	++	++	++	
Toluene	++	+	+	
Trichloroethane	++	+	+	
Trichloroethylene	++	+	+	
Xylene	++	+	+	
Acids				
Acetic acid, 25%	++	++	++	
Acetic acid, 80%	+	+	+	
Hydrofluoric acid, 25%	+	+	+	
Hydrofluoric acid, 50%	+	+	+	
Perchloric acid, 25%	-	+	-	
Phosphoric acid, up to 10%	-	+	-	
Phosphoric acid, 86%	-	+	-	
Nitric acid, 30%	-	+	-	
Nitric acid, conc.	-	-	-	
Hydrochloric acid, 15%	-	+	-	
Hydrochloric acid, 20%	-	+	-	
Sulfuric acid, 25%	+	++	+	
Sulfuric acid, 98%	-	-	-	
Trichloroacetic acid, 25%	++	+	+	
Bases				
Ammonia, 1N	+	++	+	
Ammonium hydroxide, 25%	+	+	+	
Potassium hydroxide, 32%	-	++	-	
Sodium hydroxide, 1N	+	++	+	
Sodium hydroxide, 32%	-	+	-	
Aqueous solutions				
Formaldehyde, 30%	+	+	+	
Sodium hypochlorite, 5%	-	+	-	
Hydrogen peroxide, 35%	-	++	-	
pH range				
pH 1 to 14	-	++	-	
pH 1 to 13	-	++	-	
pH 3 to 14	+	++	+	
pH 3 to 12	++	++	++	
pH 4 to 8	++	++	++	

Contact time: 24 hours at 20°C. Chemical compatibilities can be influenced by various factors. Therefore, we recommend that you confirm compatibility with the liquid you want to filter by performing a trial filtration run before you start your actual filtration.

Legend

High compatibility	++
Limited compatibility	+
Not compatible	-

