

	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	++	+	+


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.

	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	++	++	++

++ High compatibility

+ Limited compatibility

- Not compatible

 For further information, visit www.sartorius.com