

























































































































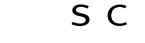









Chemical resistance of plastics





This table summarizes the resistance of plastics in Thermo Scientific™ pipettes and pipette tips to some of the most common chemicals found in laboratories. The chemical resistance table is a general guide only. Because so many factors can affect the chemical resistance of a given product, you should test the product under your specific conditions.

PP	Polypropylene
PE	Polyethylene
PVDF	Polyvinylidene fluoride
EPDM	Ethylene propylenediene monomer
FVMQ	Fluorosilicone
PEI	Polyetherimide

	Resistant (no effect)
	Limited resistance (brief exposure only)
	Not resistant
	No data available





Chemical	PP	PE	PVDF	EPDM	FVMQ	PEI
2-Chloroethanol						
Acetaldehyde						
Acetic acid, 25%						
Acetic acid, 30%						
Acetic acid, 50%						
Acetic acid, 60%						
Acetic anhydride						
Acetone						
Allyl alcohol						
Aluminum chloride						
Aluminum fluoride						
Aluminum hydroxide						
Ammonia						
Ammonia (liquid)						
Ammonium carbonate						
Amyl alcohol						
Aniline						
Barium chloride						
Benzene						
Boric acid						
Calcium chloride						

PP Polypropylene
PE Polyethylene
PVDF Polyvinylidene fluoride
EPDM Ethylene propylenediene monomer
FVMQ Fluorosilicone
PEI Polyetherimide

 Resistant (no effect)
 Limited resistance (brief exposure only)
 Not resistant
 No data available

Chemical	PP	PE	PVDF	EPDM	FVMQ	PEI
Calcium hydroxide						
Calcium sulfate						
Carbon tetrachloride						
Chlorobenzene						
Chloroform						
Chlorosulphuric acid						
Copper (II) chloride						
Dibutyl phthalate						
Dichloroethane						
Diethyl ether						
Dimethyl formamide						
Dioxane						
DMSO						
Ethanol						
Ethylene glycol						
Formaldehyde, 37%						
Formic acid (dilute)						
Furfuryl alcohol						
Glycerol						
Heptane						
Hexane						
Hydrochloric acid, 25%						
Hydrofluoric acid, 20%						
Hydrofluoric acid, 35%						
Hydrogen peroxide, 30%						
Iron (II) chloride						
Iron (III) nitrate						
Iron (III) sulfate						
Isobutanol						
Isopropanol						
Lactic acid						
Lithium bromide						
Magnesium chloride						
Magnesium nitrate						
Maleic acid						
Mercury (II) chloride						
Methanol						
Methyl ethyl ketone						
Nickel nitrate						
Nitric acid, 70%						

PP	Polypropylene
PE	Polyethylene
PVDF	Polyvinylidene fluoride
EPDM	Ethylene propylenediene monomer
FVMQ	Fluorosilicone
PEI	Polyetherimide

	Resistant (no effect)
	Limited resistance (brief exposure only)
	Not resistant
	No data available

Chemical	PP	PE	PVDF	EPDM	FVMQ	PEI
Palmitic acid						
Perchloric acid						
Phenol						
Phosphoric acid, 10%						
Polyethylene glycol						
Potassium carbonate						
Potassium chlorate						
Potassium hydroxide, 10%						
Propylene oxide						
Pyridine						
Salicylic acid						
Silver nitrate						
Sodium carbonate						
Sodium fluoride						
Sodium hydroxide, 10%						
Sodium hypochlorite, 5%						
Sulfuric acid, 50%						
Sulfuric acid, 98%						
Tannic acid						
Tetrahydrofuran						
Tin (II) chloride						
Tin (IV) chloride						
Toluene						
Trichloroacetic acid						
Triethanolamine						
Urea						
Zinc chloride						
Zinc sulfate						

Find out more at thermofisher.com/pipettes