








# Resin quick reference chart

Resin*	FEP	PPCO	PC	HDPE	PP	PS	PSF
Autoclavable	Yes	Yes	Yes	No	Yes	No	Yes
Microwavable	Marginal**	Marginal	Yes	No	Yes	No	Yes
Dry heat (oven)	Yes	No	Yes	No	No	No	Yes
Radiation	No	No	Yes	Yes	No	Yes	Yes
Freezing	Yes	Yes	Yes	Yes	No	No	Yes
Flexibility	Excellent	Moderate	Rigid	Rigid	Rigid	Rigid	Rigid
Clarity	Translucent	Translucent	Clear	Translucent	Translucent	Clear	Clear
Recycling	 OTHER	 PP	 OTHER	 HDPE	 PP	 PS	 OTHER
Noncytotoxic	Yes	Yes	Yes	Yes	Yes	Yes	Yes
Food and beverage use	Yes	Yes	Yes	Yes†	Yes	Yes	Yes

\* FEP: fluorinated ethylene propylene, PPCO: polypropylene copolymer, PC: polycarbonate, HDPE: high-density polyethylene, PP: polypropylene, PS: polystyrene, PSF: polysulfone.

\*\* The resin will absorb and retain significant amounts of heat, resulting in an unexpectedly hot surface.

† Acceptable for: nonacidic aqueous products; may contain salt, sugar, or both (pH about 5.0); dairy products and modifications; oil-in-water emulsions, high- or low-fat; moist bakery products with surface containing no free fat or oil; dry solids with surface containing no free fat or oil.

Find out more at [thermofisher.com/rigidcontainment](https://thermofisher.com/rigidcontainment)

**ThermoFisher**  
SCIENTIFIC

For Research Use or Further Manufacturing. Not for diagnostic use or direct administration into humans or animals.

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. COL05541 1217