

Chemical Resistance for Thermo Scientific Nunc Cell Culture Thermanox Coverslips

Thermo Scientific Nunc Thermanox (TMX) Coverslips are made of a polymer that is highly resistant to most chemicals. Nunc™ Thermanox™ is resistant to alcohols, aldehydes, hydro carbons, dilute acids (<10%) and dilute alkalis (<2%). Thermanox has limited resistance to chlorinated hydrocarbons, however, it is not resistant to concentrated acids or bases. It is a flexible, transparent polymer that can be sectioned using a microtome and is able to withstand high temperatures (-70°C to +150°C).

Thermanox Coverslips are available in two formats.

Rectangular:
10.5 x 22 mm
22 x 60 mm
24 x 30 mm

Round:
13 mm diameter
15 mm diameter
22 mm diameter
25 mm diameter

Thermanox Coverslips are culture treated on one side for enhanced cell attachment and growth. The treated side is packaged face up toward the label. The following tables list the chemical resistance for Thermanox Coverslips.

Chemicals having no effect	
1,2-dichloroethane	formamide
1,4-dioxane	glacial acetic acid
1-bromonaphthalene	glycerol
acetone	heptane
acetonitrile	hexyl alcohol
ammonium hydroxide (2%)	hydrochloric acid (10%)
benzene	isobutyl alcohol
butyl cellosolve	isopropanol
carbon tetrachloride	isopropyl acetate
cellosolve	isopropyl alcohol
cellosolve acetate	methanol
chloroform	methyl alcohol
cyclohexane	methyl ethyl ketone
cyclohexanol	methyl isoamyl ketone
cyclohexanone	methyl isobutyl ketone
diacetone alcohol	n-heptane
diethylene glycol	nitric acid (10%)
diethylenetriamine	n-butyl alcohol
dimethylsulfoxide	n-propyl alcohol
ethanol	sec-butyl alcohol
ethanolamine	sodium hydroxide (2%)
ethyl acetate	sulfuric acid (20%)
ethyl alcohol	tetrahydrofuran
ethylene glycol	toluene
ethylene glycol	trichloroethylene
monomethyl ether acetate	xylene

Chemicals that attack Thermanox plastic

1,1,2,2-tetrachloroethane	methyl ethyl ketone
acetic acid	methyl n-amyl ketone
acetic anhydride	methylene chloride
acetone	m-cresol
ammonium hydroxide (10%)	nitric acid (35%)
benzene	n-butyl acetate
carbon tetrachloride	n-butylamine
chloroform	n-propyl acetate
dichloroacetic acid	n-propylamine
dimethylformamide	o-chlorophenol
ethyl acetate	o-dichlorobenzene
ethylenediamine	phenol/tetrachloroethane
hexafl uoroisopropanol	sodium hydroxide (10%)
hydrochloric acid (conc.)	sulfuric acid (50%)
methyl cellosolve	tetrahydrofuran
methyl cellosolve acetate	toluene

Austria

+43 1 801 40 0

Belgium

+32 53 73 42 41

China

+86 21 68654588

Denmark

+45 4631 2000

France

+33 2 2803 2180

Germany

+49 6184 90 6940

India

+91 22 6716 2200

Italy

+39 02 02 95059 or
434-254-375

Japan

+81 3 3816 3355

Netherlands

+31 76 571 4440

Nordic/Baltic countries

+358 9 329 100

North America

+1 585-586-8800

Russia/CIS

+7 (812) 703 42 15

Spain/Portugal

+34 93 223 09 18

South America

+1 585 899 7298

Switzerland

+41 44 454 12 12

UK/Ireland

+44 870 609 9203

Other Asian countries

+852 2885 4613

Countries not listed

+49 6184 90 6940 or
+33 2 2803 2180

www.thermoscientific.com

© 2010 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

TILSPNUNCTN26 0510