

PRODUCT INFORMATION



Residual Solvents Mixture 35 (CRM)

Item No. 9004781

Supplied as: A solution in methanol

Fill volume: >1 ml

Storage: -20°C

Stability: Refer to the Certificate of Analysis

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Residual Solvents Mixture 35 (CRM) (Item No. 9004781) is a certified reference material mixture containing organic solvents. The mixture contains benzene, naphthalene, 2-propanol, n-pentane, 1,1-dichloroethene, and dichloromethane, among others. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use in method development and as a reference standard for these substances during routine analytical checks using HPLC-UV and LC-MS instrumentation. This product is intended for research and forensic applications.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/12/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



This mixture contains 35 solvents. Included are:

Item No. 9004781	Residual Solvents Mixture 35 (CRM)	
Item Name	CAS Number	Concentration (µg/ml)
Ethylene Glycol	107-21-1	5000.0
Formamide, N,N-dimethyl-	68-12-2	1500.0
Isopropyl Alcohol	67-63-0	1000.0
Pentane	109-66-0	1000.0
Ethene, 1,1-dichloro-	75-35-4	1000.0
Methylene chloride	75-09-2	1000.0
Carbon disulfide	75-15-0	1000.0
tert-Butyl methyl ether	1634-04-4	1000.0
n-Hexane	110-54-3	1000.0
Trichloromethane	67-66-3	1000.0
Ethanol, 2-methoxy-	109-86-4	1000.0
Ethane, 1,1,1-trichloro-	71-55-6	1000.0
Carbon Tetrachloride	56-23-5	1000.0
2-Propanol, 1-methoxy-	107-98-2	1000.0
Trichloroethylene	79-01-6	1000.0
1,4-Dioxane	123-91-1	1000.0
Ethanol, 2-ethoxy-	110-80-5	1000.0
Toluene	108-88-3	1000.0
Tetrachloroethylene	127-18-4	1000.0
4-Vinyl-1-cyclohexene	100-40-3	1000.0
Benzene, chloro-	108-90-7	1000.0
Ethylbenzene	100-41-4	1000.0
Styrene	100-42-5	1000.0
2-Ethoxyethyl acetate	111-15-9	1000.0
Phenol	108-95-2	1000.0
Benzene, 1,4-dichloro-	106-46-7	1000.0
Isophorone	78-59-1	1000.0
4-Phenyl-1-cyclohexene	4994-16-5	1000.0
Hexadecane	544-76-3	1000.0
Heptadecane	629-78-7	1000.0
m-Xylene	108-38-3	1000.0
o-Xylene	95-47-6	1000.0
p-Xylene	106-42-3	1000.0
Benzene	71-43-2	250.0
Naphthalene	91-20-3	250.0

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM