

PRODUCT INFORMATION



Butanes Mixture Residual Solvent (CRM)

Item No. 40824

Supplied as: A solution in 1,2,4-trimethylbenzene

Fill volume: >1 ml

Storage: -20°C

Stability: Refer to the Certificate of Analysis

Special conditions: Vortex or sonicate before use

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

Description and Data

Butanes Mixture Residual Solvent (CRM) (Item No. 40824) is a certified reference material mixture containing solvents that have been used in the extraction of cannabinoids from *Cannabis*, identified as residual solvents in *Cannabis* preparations, or included in routine threshold limits testing.¹ The mixture contains *n*-butane and isobutane. 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 threshold limits testing and in GC headspace/GC-MS applications. This product is intended for research and forensic applications.

This mixture contains 2 butanes used in the extraction of cannabinoids from *Cannabis*. Included are:

Item No. 40824	Butanes Mixture Residual Solvent (CRM)	
Item Name	CAS Number	Concentration (mg/ml)
<i>n</i> -Butane	106-97-8	2.0
Isobutane	75-28-5	2.0

Reference

1. Cannabis Laboratory Fundamentals. Opie, S.R., editor, 1st ed., Springer Cham, 325 (2021).

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, 04/05/2024

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