

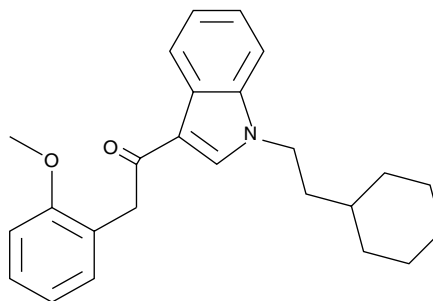
Product Information



RCS-8 (exempt preparation)

Item No. 15663

CAS Registry No.: 1345970-42-4
Formal Name: 1-(1-(2-cyclohexylethyl)-1H-indol-3-yl)-2-(2-methoxyphenyl)ethanone
Synonyms: BTM-8, SR-18
MF: C₂₅H₂₉NO₂
FW: 375.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in methanol
UV/Vis.: λ_{max}: 213, 246, 304 nm



Laboratory Procedures

For long term storage, we suggest that RCS-8 (exempt preparation) be stored as supplied at -20°C. It should be stable for at least one year.

RCS-8 (exempt preparation) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of RCS-8 (exempt preparation) in these solvents is approximately 3, 10, and 30 mg/ml, respectively.

RCS-8 (exempt preparation) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the DMF solution of RCS-8 (exempt preparation) should be diluted with the aqueous buffer of choice. RCS-8 (exempt preparation) has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

RCS-8 is a synthetic cannabinoid that has been identified as a component of several different herbal incense products. As an analog of JWH 250 (Item No. 13634), a cyclohexylethyl ring at the indole nitrogen replaces the pentyl chain. Though speculated to be less potent than JWH 250, the biological activity of RCS-8 has not been reported.

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/15663

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/26/2014