

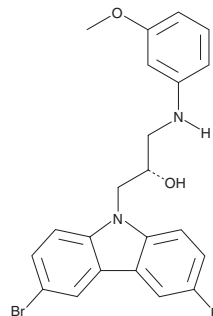
PRODUCT INFORMATION



(R)-P7C3-Ome

Item No. 34412

CAS Registry No.: 1235481-43-2
Formal Name: (αS)-3,6-dibromo-α-[[[(3-methoxyphenyl)amino]methyl]-9H-carbazole-9-ethanol
MF: C₂₂H₂₀Br₂N₂O₂
FW: 504.2
Purity: ≥98%
UV/Vis.: λ_{max}: 236, 242, 274 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

(R)-P7C3-Ome is supplied as a solid. A stock solution may be made by dissolving the (R)-P7C3-Ome in the solvent of choice, which should be purged with an inert gas. (R)-P7C3-Ome is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (R)-P7C3-Ome in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

(R)-P7C3-Ome is a hydroxylated analog of the proneurogenic and neuroprotective compound P7C3-A20 (Item No. 24201).

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, 01/03/2022

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