

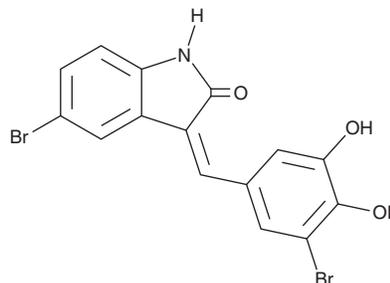
PRODUCT INFORMATION



LC3-mHTT-IN-AN1

Item No. 34306

CAS Registry No.: 486443-73-6
Formal Name: 5-bromo-3-[(3-bromo-4,5-dihydroxyphenyl)methylene]-1,3-dihydro-2H-indol-2-one
MF: C₁₅H₉Br₂NO₃
FW: 411.0
Purity: ≥95% (mixture of *cis* and *trans* isomers)
UV/Vis.: λ_{max}: 260, 372 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

LC3-mHTT-IN-AN1 is supplied as a solid. A stock solution may be made by dissolving the LC3-mHTT-IN-AN1 in the solvent of choice, which should be purged with an inert gas. LC3-mHTT-IN-AN1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LC3-mHTT-IN-AN1 in these solvents is approximately 30 mg/ml.

LC3-mHTT-IN-AN1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LC3-mHTT-IN-AN1 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. LC3-mHTT-IN-AN1 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LC3-mHTT-IN-AN1 is a linker molecule that interacts with the autophagosome protein microtubule-associated protein 1 light chain 3β (LC3B) and mutant huntingtin (Hdh).¹ It decreases Hdh^{Q7/Q140}, but not wild-type huntingtin, levels in primary mouse cortical neurons when used at a concentration of 300 nM. LC3-mHTT-IN-AN1 increases survival of *Drosophila* expressing human Hdh^{Q128}.

Reference

1. Li, Z., Wang, C., Wang, Z., *et al.* Allele-selective lowering of mutant HTT protein by HTT-LC3 linker compounds. *Nature* **575**(7781), 203-209 (2019).

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, 11/11/2022

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