

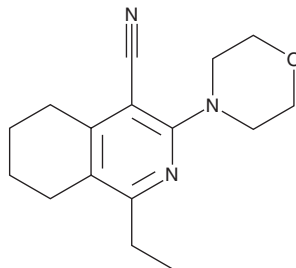
PRODUCT INFORMATION



mGluR3 Modulator-1

Item No. 45616

CAS Registry No.: 374548-18-2
Formal Name: 1-ethyl-5,6,7,8-tetrahydro-3-(4-morpholinyl)-4-isoquinolinecarbonitrile
MF: $C_{16}H_{21}N_3O$
FW: 271.4
Purity: $\geq 95\%$
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

mGluR3 modulator-1 is supplied as a solid. A stock solution may be made by dissolving the mGluR3 modulator-1 in the solvent of choice, which should be purged with an inert gas. mGluR3 modulator-1 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

mGluR3 modulator-1 is a positive allosteric modulator (PAM) of metabotropic glutamate receptor 3 (mGluR3).¹ It enhances glutamate-induced calcium mobilization in HEK293T cells expressing mGluR3 and $G_{q/15}$ ($EC_{50} = 1-10 \mu\text{M}$).

Reference

1. Potnick, J. Positive allosteric modulators of MGLUR3. **WO2014/117919A1** (2014).

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, 06/01/2026

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