

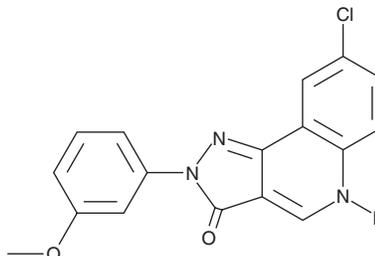
PRODUCT INFORMATION



LAU-159

Item No. 26336

CAS Registry No.: 2055050-87-6
Formal Name: 8-chloro-2,5-dihydro-2-(3-methoxyphenyl)-3H-pyrazolo[4,3-c]quinolin-3-one
MF: C₁₇H₁₂ClN₃O₂
FW: 325.8
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 290 nm
Supplied as: A crystalline 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

LAU-159 is supplied as a crystalline solid. A stock solution may be made by dissolving the LAU-159 in the solvent of choice, which should be purged with an inert gas. LAU-159 is soluble in the organic solvent DMSO and is slightly soluble in ethanol.

Description

LAU-159 is a positive allosteric modulator (PAM) of $\alpha_6\beta_3\gamma_2$ subunit-containing GABA_A receptors.¹ It selectively increases GABA-induced currents in *X. laevis* oocytes expressing $\alpha_6\beta_3\gamma_2$ over $\alpha_1\beta_3\gamma_2$, $\alpha_2\beta_3\gamma_2$, $\alpha_3\beta_3\gamma_2$, $\alpha_4\beta_3\gamma_2$, and $\alpha_5\beta_3\gamma_2$ subunit-containing receptors at 10 μ M.

Reference

1. Treven, M., Siebert, D.C.B., Holzinger, R., *et al.* Towards functional selectivity for $\alpha_6\beta_3\gamma_2$ GABA_A receptors: a series of novel pyrazoloquinolinones. *Br. J. Pharmacol.* **175**(3), 419-428 (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, 05/30/2023

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