

PRODUCT INFORMATION



ELN441958

Item No. 43214

CAS Registry No.: 913064-47-8
Formal Name: 7-chloro-2,3-dihydro-2-[3-[[9-(4-pyridinyl)-3,9-diazaspiro[5.5]undec-3-yl]carbonyl]phenyl]-1H-isoindol-1-one

MF: C₂₉H₂₉ClN₄O₂

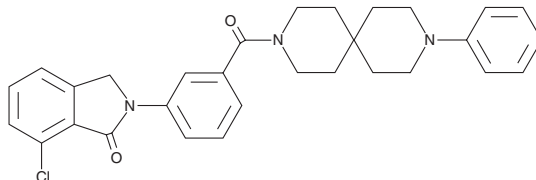
FW: 501.0

Purity: ≥98%

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

ELN441958 is supplied as a solid. A stock solution may be made by dissolving the ELN441958 in the solvent of choice, which should be purged with an inert gas. ELN441958 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

ELN441958 is a bradykinin B₁ receptor antagonist (K_i = 0.26 nM).¹ It is selective for the bradykinin B₁ receptor over a panel of greater than 100 enzymes and receptors at 10 μM but does bind to the μ-, δ-, and κ-opioid receptors and M₁ muscarinic receptor (K_is= 0.13, 0.69, 1.5, and 0.37 μM, respectively). ELN441958 inhibits calcium mobilization induced by the bradykinin B₁ receptor agonist (Des-Arg¹⁰)-kallidin (DAKD) in IMR-90 cells expressing the human bradykinin B₁ receptor (K_B = 0.12 nM). It increases tail withdrawal latencies in a rhesus monkey model of carrageenan-induced hyperalgesia when administered at doses of 3 or 10 mg/kg.

Reference

1. Hawkinson, J.E., Szoke, B.G., Garofalo, A.W., *et al.* Pharmacological, pharmacokinetic, and primate analgesic efficacy profile of the novel bradykinin B₁ receptor antagonist ELN441958. *J. Pharmacol. Exp. Ther.* **322**(2), 619-630 (2007).

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, 01/12/2026

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