

PRODUCT INFORMATION



LN6023 (hydrochloride)

Item No. 40287

Formal Name: N-(tert-butyl)-1-(3-((7-ethyl-5-oxo-5H-[1,3,4]thiadiazolo[3,2-a]pyrimidin-2-yl)amino)propyl)piperidine-4-carboxamide, dihydrochloride

MF: C₂₀H₃₂N₆O₂S • 2HCl

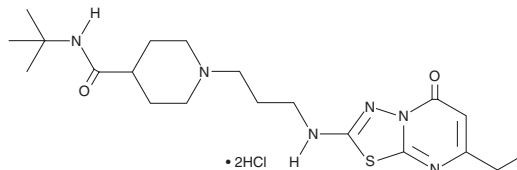
FW: 493.5

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

LN6023 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the LN6023 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. LN6023 (hydrochloride) is slightly soluble in acetonitrile. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

LN6023 is an agonist of chemokine (C-X-C motif) receptor 7 (CXCR7).¹ It induces β -arrestin recruitment in a reporter assay using HEK293T cells expressing human CXCR7 ($EC_{50} = 3.5 \mu\text{M}$) but not in HEK293T cells expressing human CXCR4. LN6023 (100 μM) reduces surface levels of P-selectin in isolated and washed human platelets.

Reference

1. Bayrak, A., Mohr, F., Kolb, K., *et al.* Discovery and development of first-in-class ACKR3/CXCR7 superagonists for platelet degranulation modulation. *J. Med. Chem.* **65**(19), 13365-13384 (2022).

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, 02/05/2024

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