

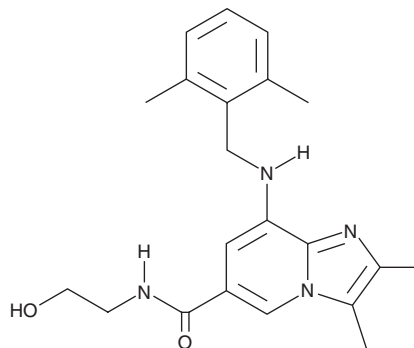
PRODUCT INFORMATION



AZD 0865

Item No. 36805

CAS Registry No.: 248919-64-4
Formal Name: 8-[[[(2,6-dimethylphenyl)methyl]amino]-N-(2-hydroxyethyl)-2,3-dimethylimidazo[1,2-a]pyridine-6-carboxamide
Synonym: Linaprazan
MF: C₂₁H₂₆N₄O₂
FW: 366.5
Purity: ≥98%
UV/Vis.: λ_{max}: 262 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

AZD 0865 is supplied as a solid. A stock solution may be made by dissolving the AZD 0865 in the solvent of choice, which should be purged with an inert gas. AZD 0865 is soluble in the organic solvent DMSO.

Description

AZD 0865 is a potassium-competitive inhibitor of the gastric H⁺/K⁺-ATPase (K_i = 46 nM).¹ It inhibits histamine- or dibutyryl-cAMP-induced acid production in isolated rabbit gastric glands (IC₅₀s = 0.28 and 0.26 μM, respectively).

Reference

1. Gedda, K., Briving, C., Svensson, K., *et al.* Mechanism of action of AZD0865, a K⁺-competitive inhibitor of gastric H⁺/K⁺-ATPase. *Biochem. Pharmacol.* **73**(2), 198-205 (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, 12/21/2022

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