

PRODUCT INFORMATION



Takeda-103A

Item No. 41027

CAS Registry No.: 865609-72-9

Formal Name: N-[(2,6-difluorophenyl)methyl]-3-[[[4-propyl-5-(4-pyrimidinyl)-4H-1,2,4-triazol-3-yl]methyl]amino]-benzamide

MF: $C_{24}H_{23}F_2N_7O$

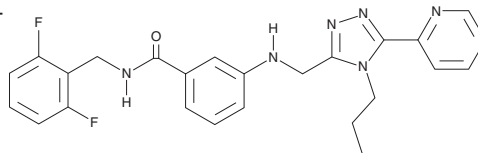
FW: 463.5

Purity: $\geq 98\%$

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

Takeda-103A is supplied as a solid. A stock solution may be made by dissolving the takeda-103A in the solvent of choice, which should be purged with an inert gas. Takeda-103A is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

Takeda-103A is an inhibitor of G protein-coupled receptor kinase 2 (GRK2).¹ It selectively inhibits phosphorylation of bovine rod outer segments by GRK2 ($IC_{50} = 54$ nM) over phosphorylation by GRK1 and GRK5 at concentrations up to 125 μM . Takeda-103A also inhibits phosphorylation of bovine rod outer segments by GRK2 containing various mutations, including GRK2^{L271M}, GRK2^{Y206S}, GRK2^{I197L}, and GRK2^{L235G} (IC_{50} s = 24, 55, 106, and 180 nM, respectively).

Reference

1. Thal, D.M., Yeow, R., Schoenau, C., *et al.* Molecular mechanism of selectivity among G protein-coupled receptor kinase 2 inhibitors. *Mol. Pharmacol.* **80**(2), 294-303 (2011).

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, 06/03/2024

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