

PRODUCT INFORMATION



TrkA Inhibitor

Item No. 45300

CAS Registry No.: 388626-12-8

Formal Name: 4-[[[(1,2-dihydro-2-oxo-3H-indol-3-ylidene)methyl]amino]benzenesulfonamide

Synonym: Tropomyosin-related Kinase A Inhibitor

MF: $C_{15}H_{13}N_3O_3S$

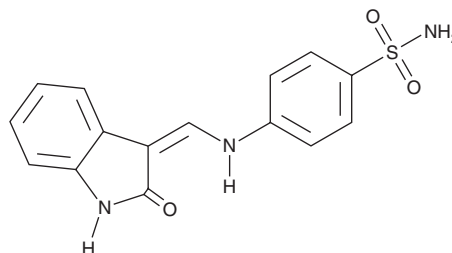
FW: 315.3

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

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

Description

TrkA inhibitor is a tropomyosin-related kinase A (TrkA) inhibitor ($IC_{50} = 0.063 \mu\text{M}$).¹ It is selective for TrkA over cyclin-dependent kinase 2 (Cdk2) and C-RAF in a cell-free assay (IC_{50} s = $>10 \mu\text{M}$ for both). TrkA Inhibitor also inhibits *T. brucei* glycogen synthase kinase 3 (GSK3) in a cell-free assay ($IC_{50} = 2.81 \mu\text{M}$).²

References

1. Wood, E. R., Kuyper, L., Petrov, K. G., *et al.* Discovery and in vitro evaluation of potent TrkA kinase inhibitors: Oxindole and aza-oxindoles. *Bioorg. Med. Chem. Lett* **14**(4), 953-7 (2004).
2. Ojo, K. K., Arakaki, T. L., Napuli, A. J., *et al.* Structure determination of glycogen synthase kinase-3 from *Leishmania major* and comparative inhibitor structure-activity relationships with *Trypanosoma brucei* GSK-3. *Mol. Biochem. Parasitol.* **176**(2), 98-108 (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, 03/12/2026

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