

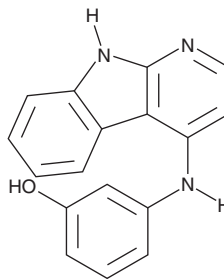
PRODUCT INFORMATION



Tilfrinib

Item No. 25647

CAS Registry No.: 1600515-49-8
Formal Name: 3-(9H-pyrido[2,3-b]indol-4-ylamino)-phenol
MF: $C_{17}H_{13}N_3O$
FW: 275.3
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

Tilfrinib is supplied as a solid. A stock solution may be made by dissolving the tilfrinib in the solvent of choice. Tilfrinib is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas. The solubility of tilfrinib in these solvents is approximately 13 and 23 mg/ml, respectively.

Description

Tilfrinib is an inhibitor of breast tumor kinase/protein tyrosine kinase 6 (BRK/PTK6; $IC_{50} = 3.15 \text{ nM}$).^{1,2} It is selective for BRK/PTK6 over HER2 ($IC_{50} = 1,300 \text{ nM}$).² Tilfrinib inhibits the growth of MCF-7, HS-578/T, and BT-549 breast cancer cells *in vitro* (GI_{50} s = 0.99, 1.02, and 1.58 μM , respectively).¹

References

1. Mahmoud, K.A., Krug, M., Wersig, T., *et al.* Discovery of 4-anilino α -carboline as novel Brk inhibitors. *Bioorg. Med. Chem. Lett.* **24**(8), 1948-1951 (2014).
2. Mahmoud, K.A., Wersig, T., Slynko, I., *et al.* Novel inhibitors of breast cancer relevant kinases Brk and HER2. *Med. Chem. Commun.* **5**, 659-664 (2014).

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, 12/19/2022

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