

PRODUCT INFORMATION



BAY-826

Item No. 41365

CAS Registry No.: 1448316-08-2

Formal Name: [3-cyano-5-[[[2,4-dimethyl-5-[6-(3-pyridinyl)-1H-imidazo[1,2-b]pyrazol-1-yl]phenyl]amino]carbonyl]phenyl] pentafluoro-sulfur

Synonyms: BAY-Tie2, BAY-Tunica Interna Endothelial Cell Kinase 2

MF: $C_{26}H_{19}F_5N_6OS$

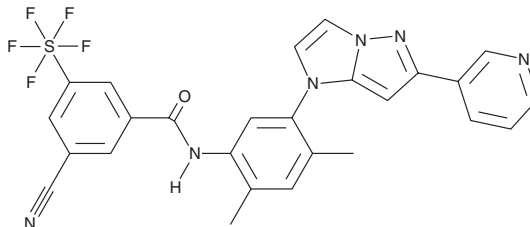
FW: 558.5

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BAY-826 is supplied as a solid. A stock solution may be made by dissolving the BAY-826 in the solvent of choice, which should be purged with an inert gas. BAY-826 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

BAY-826 is an inhibitor of tunica interna endothelial cell kinase 2 (Tie2; $IC_{50} = 0.5$ nM).¹ It inhibits Tie2 autophosphorylation in human umbilical vein endothelial cells (HUVECs; $EC_{50} = 1.3$ nM).² BAY-826 (0.1 or 1 μ M), in combination with radiation, reduces colony formation in a colony formation assay using SMA-497, SMA-540, SMA-560, and GL261 mouse glioma cells. *In vivo*, BAY 826 (100 mg/kg per day), in combination with radiation, decreases tumor volume and vascularization, as well as increases the number of CD45⁺ T cell tumor infiltrates and animal survival in an SMA-560 murine glioma model.

References

1. Sübmeier, F., Lobell, M., Grünewald, S., *et al.* Substituted phenylimidazopyrazoles and their use. *Bayer Intellectual Property*. **US2013/0190290A1** (2013).
2. Schneider, H., Szabo, E., Machado, R.A.C., *et al.* Novel TIE-2 inhibitor BAY-826 displays *in vivo* efficacy in experimental syngeneic murine glioma models. *J. Neurochem.* **140**(1), 170-182 (2016).

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, 07/17/2024

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