

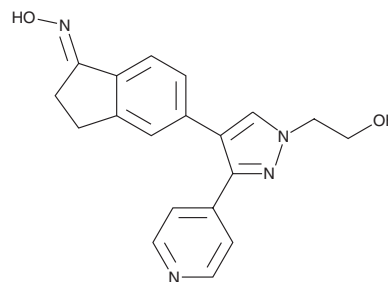
PRODUCT INFORMATION



GDC-0879

Item No. 16993

CAS Registry No.: 905281-76-7
Formal Name: 2,3-dihydro-5-[1-(2-hydroxyethyl)-3-(4-pyridinyl)-1H-pyrazol-4-yl]-1H-inden-1-one, oxime
MF: C₁₉H₁₈N₄O₂
FW: 334.4
Purity: ≥95%
UV/Vis.: λ_{max}: 267, 298 nm
Supplied as: A crystalline 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

GDC-0879 is supplied as a crystalline solid. A stock solution may be made by dissolving the GDC-0879 in the solvent of choice, which should be purged with an inert gas. GDC-0879 is slightly soluble in ethanol, DMSO, and dimethyl formamide.

Description

GDC-0879 is an inhibitor of B-Raf^{V600E} (IC₅₀ = 0.13 nM for human recombinant enzyme).¹ It is selective for B-Raf^{V600E} over a panel of 140 kinases at a concentration of 1 μM. GDC-0879 inhibits phosphorylation of ERK in MALME-3M cells (IC₅₀ = 63 nM) as well as phosphorylation of MEK1 in COLO 205 and A375 cells that express B-Raf^{V600E} (IC₅₀s = 29 and 59 nM, respectively). *In vivo*, GDC-0879 (15-200 mg/kg) inhibits tumor growth in a dose-dependent manner in an A375 mouse xenograft model.

Reference

1. Wong, H., Belvin, M., Herter, S., *et al.* Pharmacodynamics of 2-[4-[[1E]-1-(hydroxyimino)-2,3-dihydro-1H-inden-5-yl]-3-(pyridine-4-yl)-1H-pyrazol-1-yl]ethan-1-ol (GDC-0879), a potent and selective B-Raf kinase inhibitor: Understanding relationships between systemic concentrations, phosphorylated mitogen-activated protein kinase kinase 1 inhibition, and efficacy. *J. Pharmacol. Exp. Ther.* **329**(1), 360-367 (2009).

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, 12/16/2022

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