

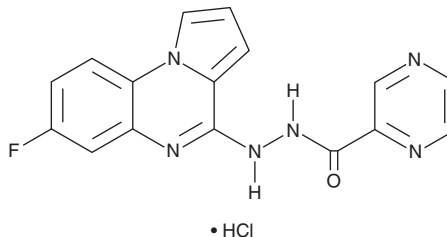
PRODUCT INFORMATION



SC144 (hydrochloride)

Item No. 29714

CAS Registry No.: 917497-70-2
Formal Name: 2-pyrazinocarboxylic acid, 2-(7-fluoropyrrolo[1,2-a]quinoxalin-4-yl)hydrazide, monohydrochloride
MF: C₁₆H₁₁FN₆O • HCl
FW: 358.8
Purity: ≥98%
UV/Vis.: λ_{max}: 269, 314 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

SC144 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SC144 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SC144 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SC144 (hydrochloride) in these solvents is approximately 5 and 2 mg/ml, respectively.

Description

SC144 is an orally bioavailable gp130 inhibitor.¹ It inhibits phosphorylation of STAT1, STAT3, and Akt induced by leukemia inhibitory factor (LIF) and IL-6, but not IFN-γ, in OVCAR-8 ovarian cancer cells when used at a concentration of 20 μM. SC144 induces cell cycle arrest at the G₀/G₁ phase in MDA-MB-435 breast and HT-29 colorectal cancer cells and inhibits proliferation of 14 cancer cell lines, including prostate, ovarian, lung, breast, and colorectal cancer cells, with IC₅₀ values ranging from 0.4 to 4 μM.² It reduces tumor growth in OVCAR-8 and MDA-MB-435 mouse xenograft models.^{1,2}

References

1. Xu, S., Grande, F., Garofalo, A., *et al.* Discovery of a novel orally active small-molecule gp130 inhibitor for the treatment of ovarian cancer. *Mol. Cancer Ther.* **12(6)**, 937-949 (2013).
2. Plasencia, C., Grande, F., Oshima, T., *et al.* Discovery of a novel quinoxalinhydrazide with a broad-spectrum anticancer activity. *Cancer Biol. Ther.* **8(5)**, 458-465 (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, 10/20/2022

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