

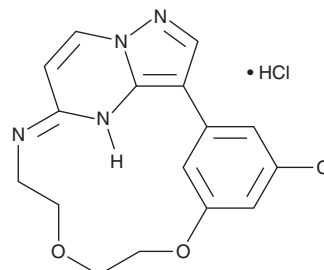
PRODUCT INFORMATION



RIPK2 Inhibitor OD36 (hydrochloride)

Item No. 42171

CAS Registry No.: 2387510-88-3
Formal Name: (1³Z)-2⁵-chloro-1⁴,1⁵-dihydro-3,6-dioxa-9-aza-1(3,5)-pyrazolo[1,5-a]pyrimidina-2(1,3)-benzenacyclononaphane, monohydrochloride
Synonyms: OD36, Receptor-interacting Serine/Threonine Kinase 2 Inhibitor OD36
MF: C₁₆H₁₅ClN₄O₂ • HCl
FW: 367.2
Purity: ≥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

RIPK2 inhibitor OD36 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the RIPK2 inhibitor OD36 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. RIPK2 inhibitor OD36 (hydrochloride) is soluble (≥10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

RIPK2 inhibitor OD36 is an inhibitor of receptor-interacting serine/threonine kinase 2 (RIPK2; IC₅₀ = 5.3 nM).¹ It also inhibits wild-type and R206H mutant activin receptor-like kinase 2 (ALK2; IC₅₀ = 47 and 22 nM, respectively).² RIPK2 inhibitor OD36 (0.5 μM) inhibits activin A-induced chondrogenic and osteogenic differentiation in ATDC-5 cells overexpressing ALK^{R206H}, an activating mutation found in patients with fibrodysplasia ossificans progressiva (FOP), which is characterized by progressive heterotopic ossification, or diffuse intrinsic pontine glioma (DIPG), a type of pediatric brain stem cancer. *In vivo*, RIPK2 inhibitor OD36 (6.25 mg/kg) reduces the number of peritoneal lavage neutrophils and lymphocytes in a mouse model of muramyl dipeptide-induced peritonitis.¹

References

1. Tigno-Aranjuez, J.T., Benderitter, P., Rombouts, F., *et al.* *In vivo* inhibition of RIPK2 kinase alleviates inflammatory disease. *J. Biol. Chem.* **289**(43), 29651-29664 (2014).
2. Sánchez-Duffhues, G., Williams, E., Benderitter, P., *et al.* Development of macrocycle kinase inhibitors for ALK2 using fibrodysplasia ossificans progressiva-derived endothelial cells. *JBMR Plus* **3**(11), e10230 (2019).

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, 12/11/2024

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