

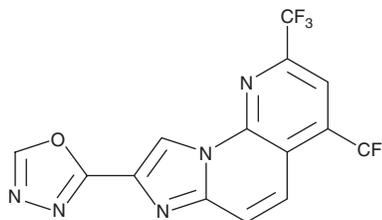
PRODUCT INFORMATION



RO8191

Item No. 20449

CAS Registry No.: 691868-88-9
Formal Name: 8-(1,3,4-oxadiazol-2-yl)-2,4-bis(trifluoromethyl)-imidazo[1,2-a][1,8]naphthyridine
Synonym: Ro 494-8191
MF: C₁₄H₅F₆N₅O
FW: 373.2
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 257, 286, 336 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

RO8191 is supplied as a crystalline solid. A stock solution may be made by dissolving the RO8191 in the solvent of choice. RO8191 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of RO8191 in these solvents is approximately 0.1 and 0.5 mg/ml, respectively.

RO8191 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RO8191 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. RO8191 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

RO8191 is an agonist of interferon (IFN) α receptor type 2 (IFNAR2) with antiviral activity (EC₅₀ = 200 nM in an anti-hepatitis C virus (HCV) replicon assay).¹ It exhibits 1:1 binding to the IFNAR2 extracellular domain, indicating RO8191 and IFNAR2 interact at the cell surface. RO8191 induces expression of IFN-stimulated genes, without induction of genes encoding inflammatory cytokines and chemokines, in HCV replicon cells, several cancer cell lines, and human primary hepatocytes. RO8191 induces expression of IFN-stimulated genes *in vivo* without induction of inflammatory cytokine and chemokine expression in mice. It also reduces HCV titer in HCV-infected humanized liver mice following oral administration.

Reference

1. Konishi, H., Okamoto, K., Ohmori, Y., *et al.* An orally available, small-molecule interferon inhibits viral replication. *Sci. Rep.* 2, 259 (2012).

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, 10/31/2022

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