

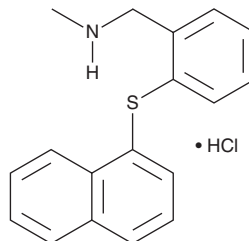
PRODUCT INFORMATION



IFN- α -IFNAR-IN-1 (hydrochloride)

Item No. 30960

CAS Registry No.: 2070014-98-9
Formal Name: N-methyl-2-(1-naphthalenylthio)-benzenemethanamine, monohydrochloride
MF: C₁₈H₁₇NS • HCl
FW: 315.9
Purity: \geq 98%
UV/Vis.: λ_{max} : 223 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

IFN- α -IFNAR-IN-1 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the IFN- α -IFNAR-IN-1 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. IFN- α -IFNAR-IN-1 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of IFN- α -IFNAR-IN-1 (hydrochloride) in these solvents is approximately 30 and 20 mg/ml, respectively.

Description

IFN- α -IFNAR-IN-1 is an inhibitor of the protein-protein interaction between the cytokine IFN- α and the IFN- α / β receptor (IFNAR; $K_d = 4 \mu\text{M}$ for binding to IFN- α).¹ It inhibits virus-induced increases in the production of IFN- α and IL-12 in bone marrow-derived plasmacytoid dendritic cells (BM-pDCs) infected with modified *Vaccinia* virus Ankara (MVA) or vesicular stomatitis virus (VSV) when used at a concentration of 18 μM . IFN- α -IFNAR-IN-1 also reduces increases in IL-12 production induced by stimulation with the toll-like receptor 9 (TLR9) agonist CpG 2216 or the TLR3 agonist poly(I:C) in the same cells.

Reference

1. Geppert, T., Bauer, S., Hiss, J.A., *et al.* Immunosuppressive small molecule discovered by structure-based virtual screening for inhibitors of protein-protein interactions. *Angew. Chem. Int. Ed. Engl.* **51**(1), 258-261 (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, 12/12/2022

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