

INCB024360 Analog

Catalog #: 27338-2

Lot #: 150203

Size: 100 mg

Structure:

CAS Registry #: 914471-09-3

Purity: ≥98%

Chemical Formula: C₉H₇ClFN₅O₂

Molecular Weight: 271.6



Description: INCB024360 analog is an orally available hydroxyamidine and inhibitor of indoleamine 2,3-dioxygenase (IDO1), with an IC₅₀ of 10 nM. It shows little activity against other related enzymes such as IDO2 or tryptophan 2,3-dioxygenase (TDO). INCB024360 analog displays potential immunomodulating and antineoplastic activities.

Synonym: 4-[(3-chloro-4-fluoroanilino)-nitrosomethylidene]-1,2,5-oxadiazol-3-amine

Appearance: White powder

Solubility: Soluble in DMSO.

Biological Activity: INCB024360 analog has been reported to restore and increase the proliferation of various immune cells including dendritic cells, NK cells, and T cells *in vitro* and *in vivo*.

Storage/Stability: Store as supplied at or below -20 °C for up to 2 years. Stable in DMSO at -80 °C for up to 6 months.

References:

1. Koblisch, HK., *et al.*, *Molec. Cancer Therapeutics*. 2010; **9**: 489-498.