

## INCB024360 Analog

**Catalog #:** 27338-1

**Lot #:** 150827

**Size:** 10 mg

**Structure:**

**CAS Registry #:** 914471-09-3

**Purity:** ≥98%

**Chemical Formula:** C<sub>9</sub>H<sub>7</sub>ClFN<sub>5</sub>O<sub>2</sub>

**Molecular Weight:** 271.6



**Description:** INCB024360 analog is an orally available hydroxyamidine and inhibitor of indoleamine 2,3-dioxygenase (IDO1), with an IC<sub>50</sub> 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.