

## Epacadostat (INCB024360)

**Catalog #:** 27339-1

**Lot #:** 151013

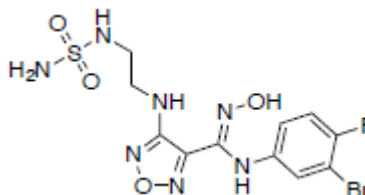
**Size:** 10 mg

**CAS Registry #:** 1204669-58-8

**Structure:**

**Purity:** ≥98%

**Chemical Formula:** C<sub>11</sub>H<sub>13</sub>BrFN<sub>7</sub>O<sub>4</sub>S



**Molecular Weight:** 438.2

**Description:** Epacadostat (INCB024360) is an orally available hydroxylamine 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). Epacadostat displays potential immunomodulating and antineoplastic activities.

**Synonym:** INCB024360, INCB24360

**Appearance:** White powder

**Solubility:** Soluble in DMSO.

**Biological Activity:** Epacadostat 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.