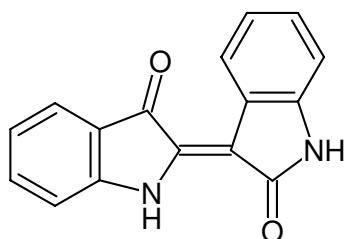


**Indirubin**

Inhibitor of Aurora Kinases
 Indigo red
 Indigopurpurin
 Indigo red

Cat. No.	Amount
NU-964	1 mg



Structural formula of Indirubin

For research use only!**Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₁₆H₁₀N₂O₂**Molecular Weight:** 262.26 g/mol**CAS#:** 906748-38-7**Purity:** ≥ 95 % (HPLC)**Form:** solid**Applications:****For research use only!****Description:**

Aurora kinases are serine/threonine kinases. They play a role in cell division by controlling chromatid segregation. Aurora kinase A is implicated in important processes during mitosis including centrosome and microtubule activities. The substrate protein Aurora A is phosphorylated at threonine 248. Phosphorylation of histone H3 on serine 10 is required for proper execution of mitosis and precedes Aurora A dependent. Indirubin and indirubin-derivatives inhibit the proliferation of a variety of cells. Thus Aurora kinases are an attractive target for anticancer therapy.

Selected References:

Muzinga Ndolo *et al.* (2017) Characterization of the indirubin derivative LDD970 as small molecule Aurora kinase A inhibitor in human colorectal cancer cells. *Immune network* **17** (2):110.

Conner (2015) Oncolytic virus and aurora kinase inhibitor for treatment of cancer. *PCT Int. Appl.* WO 2015173544 A1 20151119

Brown (2013) Compositions and methods to improve the therapeutic benefit of indirubin and analogs, including mesoindigo. *PCT Int. Appl.* WO 2013142817 A2 20130926.

Mahindra *et al.* (2010) Novel therapeutic targets for multiple myeloma. *Future in Oncology* **6** (3):407.

Myrianthopoulos *et al.* (2007) An integrated computational approach to the phenomenon of potent and selective inhibition of Aurora kinases B and C by a series of 7-substituted indirubins. *J. Med. Chem.* **50**:4027.

Bolanus-Garcia (2005) Aurora kinases. *Int. J. Biochem. & Cell Biol.* **37**:1572.