

LAQ824 (NVP-LAQ824, Dacinostat)

Catalog #: 27760-1

Lot #: 230208

Size: 50 mg

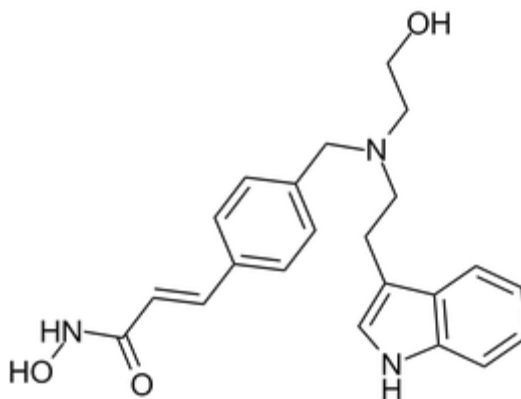
Structure:

Cas Registry: 404951-53-7

Purity: ≥98%

Chemical Formula: C₂₂H₂₅N₃O₃

Molecular Weight: 379.46 Da



Background: A recent study shows that LAQ824 induces chromatin changes at the level of the IL-10 gene promoter that lead to enhanced recruitment of the transcriptional repressors HDAC11 and PU.1 and inhibits IL-10 production in BALB/c murine macrophages.

Appearance: A solid

Description: LAQ824 (also known as NVP-LAQ824 or Dacinostat), a derivative of 4-aminomethylcinnamic hydroxamic acid, is a novel and potent inhibitor of histone deacetylase (HDAC) that inhibits the activity of HDAC with 50% inhibition concentration IC₅₀ value of 0.03 μM. LAQ824 has been found to inhibit the growth of a variety of cancer cell lines, including colon cancer H1299 and HCT116 cells, breast cancer MDA435 cells, prostate cancer DU145 and PC3 cells and non-small cell lung cancer A549 cells, with IC₅₀ value < 1 μM and induce apoptosis in human breast cancer SKBR-3, BT-474 and MB-468 cells. LAQ824 also dose- and time-dependently inhibits the growth of multiple myeloma cells.

Solubility: insoluble in H₂O; insoluble in EtOH; ≥17.45 mg/mL in DMSO

Storage/Stability: Store at or below -20°C. Solid form is stable at least 12 months from date of receipt, when stored as directed. Do not store aqueous solutions for more than one day.

Reference:

1. Catley L, et al. Blood. 2003 Oct 1;102(7):2615-22.