

Product Information

Description

Trichostatin A (TSA) is a potent inhibitor of histone deacetylase (HDAC) as well as an antifungal antibiotic with cytostatic and differentiating properties that noncompetitively and reversibly inhibits HDAC, at low nanomolar concentrations, in both cultured mammalian cells and fractionated cell nuclear extracts.

Target:

DNA Damage/DNA Repair

Pathway:

HDAC

Chemical Name:

(2E,4E,6R)-7-[4-(dimethylamino)phenyl]-N-hydroxy-4,6-dimethyl-7-oxohepta-2,4-dienamide

Molecular Formula:

C₁₇H₂₂N₂O₃

Supplied As:

Solid

Solubility:

insoluble in H₂O; ≥15.12 mg/mL in DMSO; ≥16.56 mg/mL in EtOH with ultrasonic. For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while.

Stability:

At least 6 months at -20°C. Avoid freeze/thaw cycles.

Storage:

-20°C

MW:

302.37 Da

Cas No.

58880-19-6

Purity:

≥98% by HPLC

Chemical Structure

