

Product Information

Description

Potent MDM2 inhibitor with an IC₅₀ = 6 nM in binding assay and 30 nM in cancer cell proliferation assay. Induces p53 stabilisation, cell cycle arrest and apoptosis in cancer cells expression wildtype p53. Displays inhibition of tumor growth in the SJSA1 tumor xenograft model. Also inhibits MDR-1 at high concentrations.

Molecular Formula:

C₃₁H₂₉Cl₂F₂N₃O₄

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

616.5 Da

Cas No.

1229705-06-9

Chemical Structure

