

Product Information

Description

Alvespimycin HCl, also known as 17-DMAG, is a potent, water-soluble HSP90 inhibitor with an IC_{50} value of 62 ± 29 nM. Alvespimycin has diverse anti-tumor actions and has potential in treating a number of different types of cancers, including ovarian, prostate, melanoma, non-small-cell lung cancer, and several leukemias (chronic lymphocytic, acute myeloid). This compound also suppresses inflammation by interfering with signaling through the NF- κ B pathway.

Target:

HSP90

Pathway

Metabolism

Chemical Name:

[(3R,5S,6R,7S,8E,10S,11S,12Z,14E)-21-[2-(dimethylamino)ethylamino]-6-hydroxy-5,11-dimethoxy-3,7,9,15-tetramethyl-16,20,22-trioxo-17-azabicyclo[16.3.1]docosa-1(21),8,12,14,18-pentaen-10-yl] carbamate;hydrochloride

Molecular Formula:

 $C_{32}H_{48}N_4O_8 \cdot HCl$

Supplied As:

Solid

Solubility:

≥ 26.2 mg/mL in DMSO; insoluble in EtOH; ≥ 3.04 mg/mL in H_2O For obtaining a higher solubility, please warm the tube at $37^\circ C$ and shake it in the ultrasonic bath for a while. Long-term storage in solution is not recommended. Not stable in diluted solutions.

Stability:

At least 6 months at $-20^\circ C$. Avoid freeze/thaw cycles.

Storage:

 $-20^\circ C$

MW:

653.21 Da

Cas No.

467214-21-7

Purity:

 $\geq 99\%$

Chemical Structure

