Necrostatin-5

Necroptosis inhibitor

Necroptosis inhibitor. In contrary to the mechanism of necrostatin-1 (Prod.No. BML-AP309) it was shown that necrostatin-5 inhibits RIP1 kinase indirectly. A potent inhibitor of immunoprecipitated, but not recombinant RIP1.

The necrostatins inhibit non-apoptotic cell death pathway by inhibition of the death domain receptor-associated adaptor kinase RIP1. There are three different mechanisms: T-loop dependent inhibition by necrostatin-1, partially T-loop independent inhibition by necrostatin-3 and indirect inhibition of RIP1 by necrostatin-5.

Citations: 3

View Online »

Ordering Information

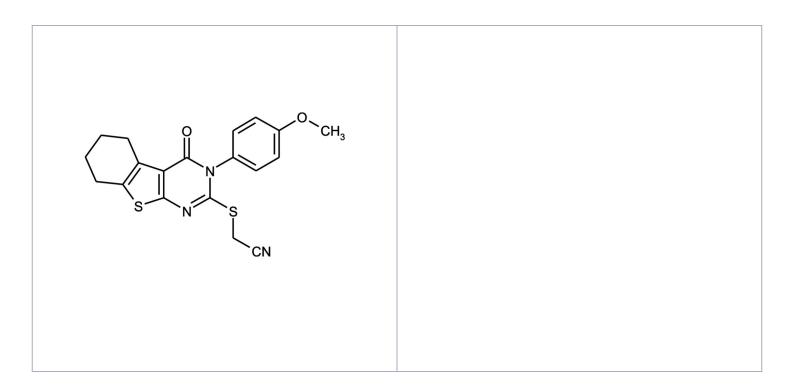
Order Online »

ALX-430-169-M025

25mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Nec-5, 3-p-Methoxyphenyl-5,6-tetramethylenothieno[2,3-

d]pyrimidin-4-one-2-mercaptoethylcyanide

Appearance White to off-white solid.

CAS 337349-54-9

Couple Target RIP

Couple Type Inhibitor

Formula $C_{19}H_{17}N_3O_2S_2$

MW 383.5

Purity ≥98% (HPLC, TLC)

Soluble in DMSO (>10mg/ml).

Last modified: May 29, 2024