

BAY-11-7082

Catalog #: 27229

Lot #: 120628

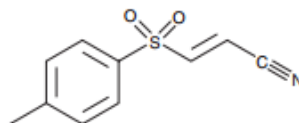
Size: 10 mg

Structure:

CAS Registry #: 19542-67-7

Purity: ≥98%

Chemical Formula: C₁₀H₉NO₂S



Molecular Weight: 207.3

Description: BAY-11-7082 is a selective and irreversible inhibitor of NF-κB activation by blocking TNF-α-induced phosphorylation of IκB-α without affecting constitutive IκB-α phosphorylation. It also inhibits the TNF-α-induced surface expression of adhesion molecules ICAM-1, VCAM-1, and E-selectin in human endothelial cells.

Appearance: A crystalline solid

Solubility: Soluble in ethanol, DMSO and DMF, purged with an inert gas. Solubility in ethanol is ~ 0.20 mg/ml, and in DMSO and DMF is 25 mg/ml. Do not store aqueous solutions for more than one day.

Biological Activity: BAY-11-7082 inhibits TNF-α-induced surface expression of adhesion molecules ICAM-1, VCAM-1, and E-selectin with IC₅₀ values of 5-10 μM.

Storage/Stability: Store at or below -20 °C for up to two years.

Quality Control: The purity was determined by HPLC analysis.

Reference(s):

1. Pahl, H., *et al.*, *Oncogene*. 1999;**18**:6853-6866.
2. Gilmore, TD, *et al.*, *Oncogene*. 1999;**18**:6842-6844.
3. Karin, M., *et al.*, *J. Biol. Chem.* 1999;**274(39)**:27339-27342.
4. Pierce, J.W., *et al.*, *J. Biol. Chem.* 1997;**272(34)**:21096-21103.