

## U0126

**Catalog #:** 27012

**Lot #:** 181219

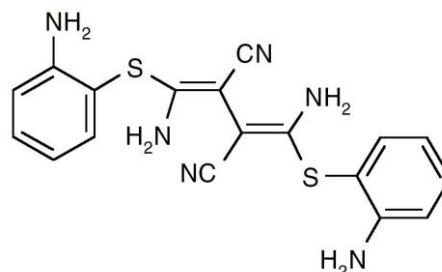
**Size:** 25 mg

**Structure:**

**CAS Registry #:** 109511-58-2

**Purity:** ≥99%

**Chemical Formula:** C<sub>18</sub>H<sub>16</sub>N<sub>6</sub>S<sub>2</sub>



**Molecular Weight:** 380.49

**Description:** U0126 is a selective inhibitor of mitogen-activated protein kinase kinases, MEK-1 and MEK-2. It is a weak inhibitor of PKC, Raf, ERK, JNK, MEKK, MKK-3, MKK-4/SEK, MKK-6, Abl, Cdk2 and Cdk4.

**Appearance:** White to off-white crystalline solid; granular or powder

**Solubility:** Soluble in DMSO at 200 mg/ml. It is poorly soluble in ethanol and water. Buffers, serum or other additives may increase or decrease the aqueous solubility.

**Biological Activity:** U0126 antagonizes AP-1 transcriptional activity via noncompetitive inhibition of the dual specificity kinase MEK with IC<sub>50</sub> of 0.07 μM for MEK1 and 0.06 μM for MEK2.

**Storage/Stability:** Store at or below -20°C.

**Quality Control:** The purity was determined by HPLC analysis.

### References:

1. Favata, M.R., *et al.*, *J. Biol. Chem.* 1998; **273**: 18623 – 18632.
2. DeSilva, D.R., *et al.*, *J. Immunol.* 1998; **160**: 4175 – 4181.
3. Scherle, P.A., *et al.*, *J. Immunol.* 1998; **161**: 5681 – 5686.
4. Duncia, J.V., *et al.*, *Bioorg Med Chem Lett.* 1998 Oct 20; **8(20)**: 2839 – 2844.



## Data Sheet