Tyrphostin 51

EGFR inhibitor

A potent inhibitor of EGF receptor kinase activity (IC $_{50}$ =0.8 μM). Inhibition is mixed competitive with ATP and substrate.

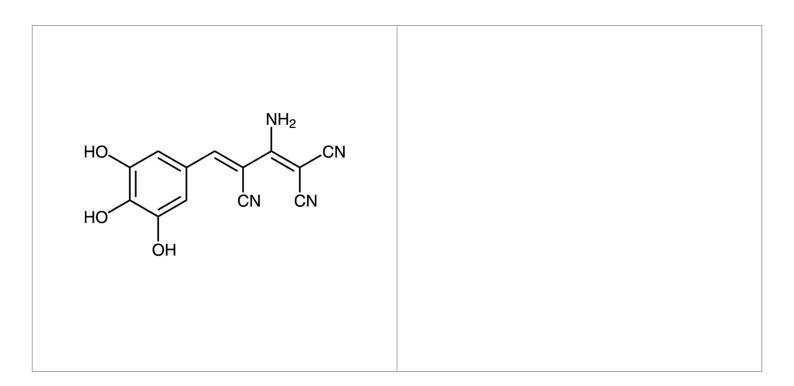
Ordering Information

Order Online »

BML-EI189-0005 5mg

Manuals, SDS & CofA

View Online »



Handling & Storage

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

at-20°C for up to 1 year. Product is subject to oxidation. Make solutions fresh daily.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Brown solid.

CAS 122520-90-5

Couple Target EGFR

Couple Type Inhibitor

Formula $C_{13}H_8N_4O_3$

MW 268.2

Purity ≥99%

Solubility Slightly soluble in water.

Last modified: May 29, 2024

