

AG-370

PDGF receptor inhibitor

Member of the tyrphostin family of tyrosine kinase inhibitors. Selective inhibitor of the PDGF receptor kinase ($IC_{50}=20\mu\text{M}$) vs. the EGF receptor kinase ($IC_{50}=820\mu\text{M}$). Inhibits PDGF-induced mitogenesis in human bone marrow fibroblasts.

Citations: 2

[View Online »](#)

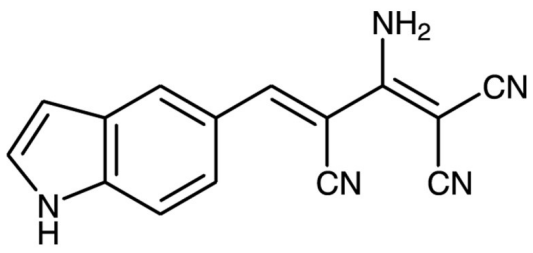
Ordering Information

[Order Online »](#)

BML-EI229-0002	2mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Store tightly sealed in the dark. Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Tyrphostin AG 370, Tyrphostin B7, 2-Amino-4-(1H-indol-5-yl)-1,1,3-tricyanobuta-1,3-diene
Appearance	Yellow solid.
CAS	134036-53-6
Couple Target	PDGF receptor
Couple Type	Inhibitor
Formula	$C_{15}H_9N_5$
MW	287.3
MeltingPoint	300-302°C
Purity	≥99% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (2mg/ml, warm).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com