

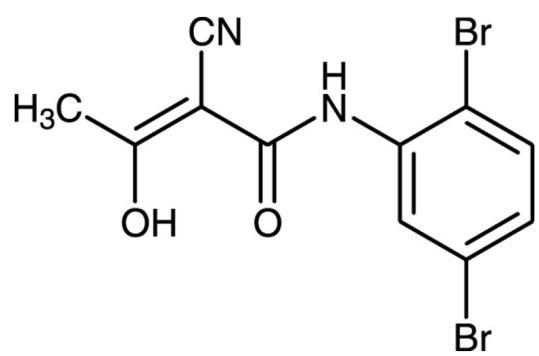
LFM-A13

BTK inhibitor

Bruton's tyrosine kinase (BTK) inhibitor $IC_{50} = 2.5 \mu M$. IC_{50} 's for JAK-1, JAK-2, JAK-3, SYK, HCK, EGFR kinase, IR kinase all $>300 \mu M$

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. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Yellow solid .
CAS	62004-35-7, 244240-24-2
Couple Target	BTK
Couple Type	Inhibitor
Formula	$C_{11}H_8Br_2N_2O_2$
MW	360.0
MeltingPoint	145-146°C
Purity	≥97% (TLC)
Solubility	Soluble in DMSO (10mg/ml) or ethanol (0.3mg/ml).

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-fr@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