

STA-21

STAT inhibitor

STA-21 is a novel Stat3-selective inhibitor. It inhibits Stat3 dimerization and DNA binding as well as Stat3-dependent luciferase reporter activity.

Citations: 4

[View Online »](#)

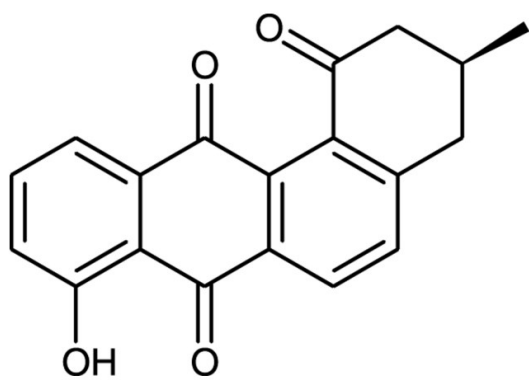
Ordering Information

[Order Online »](#)

BML-EI351-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(S)-Ochromycinone, Deoxytetrangomycin
Appearance	White to yellow solid.
CAS	28882-53-3
Couple Target	STAT
Couple Type	Inhibitor
Formula	$C_{19}H_{14}O_4$
MW	306.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (20mg/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