

Auranofin

Anti-arthritic agent

Gold derivative. Widely used as an anti-arthritic drug. Inhibits various leukocyte activation pathways at multiple sites. Inhibits release of inflammatory mediators from human basophils, pulmonary mast cells and macrophages. Also inhibits human neutrophil 5-lipoxygenase. Inhibits I κ B kinase (IKK) by modifying Cys¹⁷⁹ of the IKK β subunit. Efficient inducer of mitochondrial membrane permeability transition pore in the presence of calcium ions, potentially referable to its inhibition of mitochondrial thioredoxin reductase.

Citations: 30

[View Online »](#)

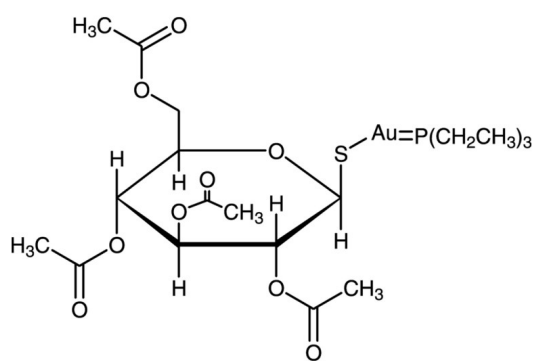
Ordering Information

[Order Online »](#)

BML-EI206-0100	100mg
----------------	-------

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, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	SKF 39162, 1-Thio-β-D-glucopyranosatotriethyl phosphine gold-2,3,4,6-tetraacetate
Appearance	White solid.
CAS	34031-32-8
Couple Target	IKK, Lipoxygenase
Couple Type	Inhibitor
Formula	$C_{20}H_{34}AuO_9PS$
MI	14: 878
MW	678.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (5mg/ml) or 100% ethanol (4mg/ml).



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