

Cafestol

GST activator

Inducer of glutathione S-transferases. Has chemoprotective activity, reducing the genotoxicity of several carcinogens.

Citations: 5

[View Online »](#)

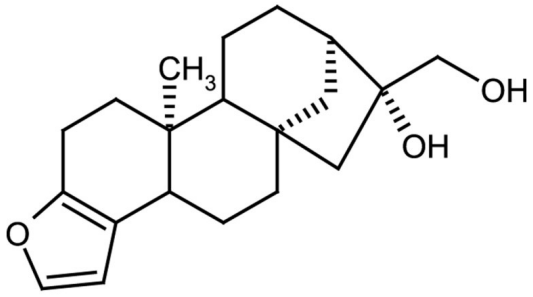
Ordering Information

[Order Online »](#)

ALX-350-220-M050	50mg
ALX-350-220-46190	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white solid.

CAS 469-83-0

Couple Target GST

Couple Type Activator

Formula $C_{20}H_{28}O_3$

MI **14:** 1634

MW 316.4

Purity ≥98% (UHPLC)

Solubility Soluble in DMSO, 100% ethanol or ethyl acetate.

Source Isolated from the unsaponifiable fraction of petroleum ether extract of coffee beans.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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