

Timosaponin AIII

mTOR inhibitor

Induces autophagy in HeLa cells followed by apoptotic cell death (IC_{50} =10 μ M). This cellular response has been found to be mediated *via* inhibition of mTORC1 and induction of ER stress (IC_{50} for BT474 cells=2.5 μ M and MDAMB231 cells=6 μ M). The pro-apoptotic response is selective for tumor cells over normal cells but autophagy is induced in both.

Citations: 3

[View Online »](#)

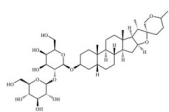
Ordering Information

[Order Online »](#)

BML-AP504-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 6 months after receipt when stored at -20°C. Solutions can be stored at -20°C for up to 3 months.
Handling	Solutions can be stored at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off white solid.
CAS	41059-79-4
Couple Target	mTOR
Couple Type	Inhibitor
Formula	$C_{39}H_{64}O_{13}$
MW	740.9
Purity	≥96% (HPLC)
Solubility	Soluble in DMSO (20 mg/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