

ONO-RS-082

Phospholipase inhibitor

Phospholipase A₂ inhibitor (IC₅₀=7 µM for purified guinea pig lung PLA₂). Inhibits epinephrine-stimulated thromboxane production in human platelets (86% at 3.5µM). Also blocks insulin secretion in pancreatic islets in response to glucose. Disrupts membrane tubule-mediated reassembly and maintenance of the Golgi complex.

Citations: 6

[View Online »](#)

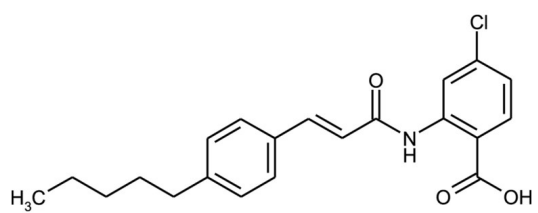
Ordering Information

[Order Online »](#)

BML-ST320-0020	20mg
BML-ST320-0100	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 2 years after receipt when stored at +20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light. Do not freeze.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-(p-Amylcinnamoyl)amino-4-chlorobenzoic acid
Appearance	Off-white to yellow solid.
CAS	99754-06-0
Couple Target	Phospholipase, PLA
Couple Type	Inhibitor
Formula	$C_{21}H_{22}ClNO_3$
MW	371.9
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml, warm) or 100% ethanol (10mg/ml, warm).

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)