

Z-Phe-Arg-pNA . hydrochloride

Cathepsin substrate

Chromogenic substrate for Cathepsins B, K, L, and S, papain, trypsin, and plasma kallikrein. This substrate is likely to be cleaved by other cathepsins (see Z-Phe-Arg-AMC). Release of pNA is monitored at 405-410 nm (pNA $\epsilon_{410}=8800 \text{ M}^{-1}\text{cm}^{-1}$). Useful for inhibitor screening and kinetic analysis.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-130-M005	5mg
------------------	-----

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

Alternative Name Z-FR-pNA . HCl

Appearance White to off-white powder.

CAS 59188-54-4

Formula $C_{29}H_{33}N_7O_6 \cdot HCl$

Formulation Lyophilized

MW 575.6 . 36.5

Purity ≥99% (TLC)

Sequence Z-Phe-Arg-pNA (pNA=p-Nitroanilide)

Solubility Soluble in methanol (50mg/ml) and DMSO.

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-
fr@enzolifesciences.com](mailto:info-fr@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)