

Bactenecin-5

precursor peptide

Bac5 is a 5-kDa proline-arginine rich antibiotic with a high degree of homology with PR39, an inhibitor of the 20S and 26S proteasome. Recently bactenecin 5 has been demonstrated to inhibit the chymotrypsin-like activity of the 20S proteasome.

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Bac5-GRR

Appearance White powder.

MW 5517.8

Purity ≥95% (HPLC)

Sequence H-Arg-Phe-Arg-Pro-Pro-Ile-Arg-Arg-Pro-Pro-Ile-Arg-Pro-Pro-Phe-Tyr-Pro-Pro-Phe-Arg- Pro-Pro-Ile-Arg-Pro-Pro-Ile-Phe-Pro-Pro-Ile-Arg-Pro-Pro-Phe-Arg-Pro-Pro-Leu-Gly-Pro-Phe-Pro-Gly-Arg-Arg-OH

Solubility Soluble in aqueous solution.

Last modified: May 29, 2024



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