

Substance P

Neuropeptide

Neuropeptide inducing Ca^{2+} mobilization. Activates G_1 -Proteins directly. Induces superoxide production in neutrophils.

Substance P (SP) was first discovered in 1931 by von Euler and Gaddum and is the best characterized member of the neurokinins. This class of peptide neurotransmitters includes neurokinin A (NK-A) and B (NK-B), and are characterized by the common C-terminal sequence Phe-X-Gly-Leu-Met- NH_2 . The receptors for SP, NK-A and NK-B, designated NK_1 , NK_2 and NK_3 respectively, are members of the G protein-coupled receptor superfamily. All three receptors have been shown to stimulate inositol phospholipid hydrolysis. SP is involved in pain, smooth muscle contraction, secretion from exocrine and endocrine glands, inflammatory response including vasodilation, plasma extravasation and mobilization of immune system cells.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-167-003-M001	1mg
ALX-167-003-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

CAS 33507-63-0

Formula $C_{63}H_{98}N_{18}O_{13}S$

Formulation Lyophilized

MI 14: 8863

MW 1347.6

Purity ≥97% (HPLC)

Sequence H-Arg-Pro-Lys-Pro-Gln-Gln-Phe-Phe-Gly-Leu-Met-NH₂

Solubility Soluble in DMSO or water.



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)