

BAM (8-22)

BAM (8-22) is a opioid peptide fragment of the bovine adrenal medulla (BAM22) neuropeptide that is involved in inflammation, pain, and morphine tolerance. Induces itch and nociceptive sensations. Lacks the met-enkephalin motif of BAM22, hence lacking affinity for opioid receptors. Selective agonist of the sensory neuron specific receptors SNSR3 and SNSR4 (EC50=28 and 14nM, respectively).

Materials Provided

Catalog Number:	92-1215
Quantity Supplied:	1 vial(s), 1 mg per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	1971.22 g/mol
Source:	Synthetic
Purity:	≥ 81%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	1 mM
Reconstitution Volume:	0.507 mL
Reconstitution Solvent:	H ₂ O

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com