

Product Name:	Caerulein	
Catalog Number:	AS-24252 (1 mg)	Lot Number: See label on vial
Sequence:	Pyr-Gln-Asp-Tyr(SO3H)-Thr-Gly-Trp-Me Pyr-QD-Y(SO3H)-TGWMDF-NH2 (1-lett	,
Molecular Weight:	1353.4	
% Peak Area by HPLC:	Area by HPLC: ≥ 95	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Reconstitute by adding 100 µl 1%NH₄OH to 1 mg Caerulein peptide. Dilute this peptide solution to approximately 1 mg/ml (or more dilute) with a buffer such as PBS or another buffer; aliquot and store at -20C.

Storage: Caerulein peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20° C or lower. Reconstituted peptide can be aliquoted and stored at -20° C or lower.

Description: Caerulein, a decapeptide analog of the potent pancreatic secretagogue cholecystokinin, stimulates gastric, biliary, and pancreatic secretion. Caerulein injections cause acute pancreatitis in mice. Ref: Wabnitz, PA. et al. *Rapid Commun. Mass Spectrom.* **13**, 2498 (1999); Loguercio, C. et al. *Digestion* **48**, 128 (1991).

Related Products:

Name		Size
[Thr28, Nle31]-Cholecystokinin (25-33), sulfated (RD-Y(SO3H)-TGW-Nle-DF-NH2)	AS-22944	1 mg
Cholecystokinin (26-33), CCK Octapeptide, sulfated (D-Y(SO3H)-MGWMDF-NH2)	AS-20741	1 mg
Cholecystokinin (26-33), CCK8 (DYMGWMDF-NH2)	AS-20739	1 mg

For Research Use Only