

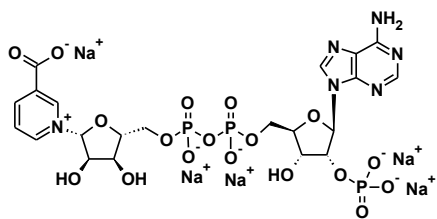
Product Information Sheet

Ordering Information

Product Number:	20999
Product Name:	NAADP [Nicotinic acid adenine dinucleotide phosphate sodium salt]
Unit Size:	1 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 855.31



Chemical Structure:

Soluble in: Water

Application Notes

NAADP is a calcium mobilizing agent. It initiates calcium release via type 1 ryanodine receptor (RyR1) activation. NAADP also regulates calcium release from intracellular stores distinct from the endoplasmic reticulum (ER), and displays affinity for two-pore channels (TPCs), which release calcium from acidic organelles.