

NAADP-AM

Catalog Number: 20997 20998, 21000,

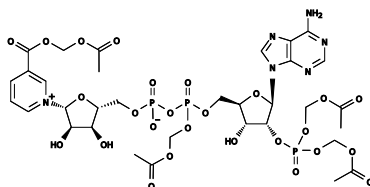
Unit Size: 2X50 ug 250 ug, 1 mg,

Product Details

Storage Conditions	Freeze (< -60 °C), Minimize light exposure,
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid light yellow
Molecular Weight	1032.64
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

NAADP (nicotinic acid adenine dinucleotide phosphate) is a recently discovered second messenger. NAADP is negatively charged (to prevent it from leaking out of cells). The negative charge of NAADP makes it well retained in cells, but imposes difficulties in loading NAADP into live cells. NAADP must be injected, loaded via liposomes, or electroporated. These invasive techniques are highly technically demanding and are possible only in certain single cell preparations. NAADP-AM has been reported as a cell-permeant analog of NAADP. NAADP-AM can be taken into cells and induces NAADP-mediated calcium signaling. NAADP-AM can be used in a wide range of systems and will greatly facilitate research into the role of NAADP in biological research. The NAADP-AM consists of three NAADP AM esters with one major component and two minor AM esters.