

NAADP-AM

Catalog Number: 20997 20998, 21000,

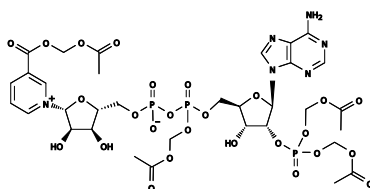
Unit Size: 2X50 ug 250 ug, 1 mg,

Product Details

| | |
|--------------------|---|
| Storage Conditions | Freeze (< -60 °C), Minimize light exposure, |
| Expiration Date | 3 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.