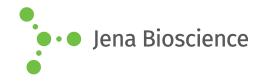
DATA SHEET





AMP-α-F

(ApF)

Adenosine-5'-(α -fluoro)-monophosphate, Sodium salt Adenosine-5'-(1-fluoro)-monophosphate, Sodium salt

Cat. No.	Amount
NU-943-5	5 mg
NU-943-25	25 mg

Structural formula of AMP-α-F

For general laboratory use.

Shipping: shipped on gel packs **Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery **Molecular Formula:** C₁₀H₁₃FN₅O₆P (free acid) **Molecular Weight:** 349.21 g/mol (free acid)

Exact Mass: 349.06 g/mol (free acid)

CAS#: 19375-33-8 (free acid), 15503-75-0 (sodium salt)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{max} 259 nm, ϵ 15.4 L $mmol^{\text{--}1}$ (Tris-HCl

pH 7.5)

Applications:

Substrate for snake venom phosphodiesterase [1,2]

Substrate for Fhit proteins [3]

Inhibitor of adenlyate kinase [4]

Selected References:

[1] Baranowski *et al.* (2016) A fluorescent HTS assay for phosphohydrolases based on nucleoside 5'-fluorophosphates: its application in screening for inhibitors of mRNA decapping scavenger and PDE-I. *Org. Biomol. Chem.* **14 (20)**:4595.

[2] Baranowski *et al.* (2015) Synthesis of fluorophosphate nucleotide analogues and their characterization as tools for ¹⁹F NMR studies. *J. Org. Chem.* **80** (8):3982.

[3] Guranowski *et al.* (2008) Fhit proteins can also recognize substrates other than dinucleoside polyphosphates. *FEBS Lett.* **582 (20)**:3152.

[4] Scoblov et al. (1996) Modified nucleotides as substrates and inhibitors of adenylate kinase from different sources. FEBS Lett. **395 (2-3)**:283.