

AMP- α -F (25 mg)

Substrate for snake venom phosphodiesterase. Substrate for Fhit proteins.
Inhibitor of adenylate kinase.

Ordering Information

[Order Online »](#)

JBS-NU-943-25	25mg
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Appearance white to off-white solid

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

Couple Type Modified nucleotides

Formula $C_{10}H_{13}FN_5O_6P$ (free acid)

MW 349.21 (free acid)

Purity $\geq 95\%$ (HPLC)

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com