

6-Fu-ATP- γ -S

NTP analogue for chemical genetics

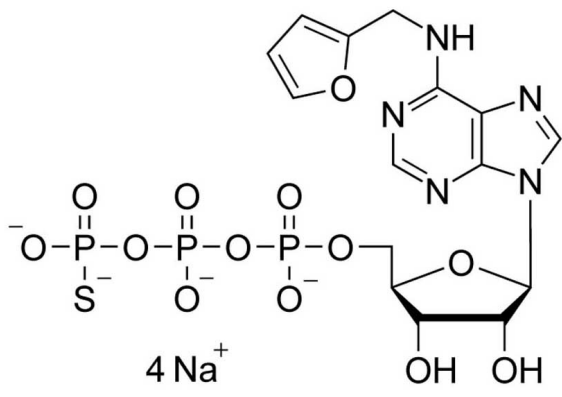
6-Fu-ATP- γ -S is useful for specific interaction with modified receptor proteins and can be used in studies on kinase-substrate relationships (chemical genetics). The related structures 6-Bn-ATP- γ -S (Prod. No. [BLG-B072](#)) and 6-PhEt-ATP- γ -S (Prod. No. [BLG-P026](#)) are offered as well.

BLG-F008-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information [Order Online »](#)

BLG-F008-05	5 μ mol
BLG-F008-25	5x5 μ mol

Manuals, SDS & CofA [View Online »](#)



Handling & Storage

Use/Stability	6-Fu-ATP- γ -S is most stable when stored as aqueous solution in the freezer (-20° Celsius necessary, -80° recommended), however, at ambient temperature the compound slowly starts to decompose. Thus, in order to maintain its original high quality it is recommended to allow thawing only before using the product. If you will not use up the vial with one application, please aliquot the contents of the vial in order to avoid repeated freeze/thaw cycles for the rest. When preparing such aliquots be sure to operate quickly and to freeze the vial again as soon as possible.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	N6-Furfuryladenosine-5'-O-(3-thiotriphosphate) . tetrasodium salt
Appearance	10 mM aqueous solution
CAS	1400539-83-4
Couple Type	Modified nucleotides
Formula	$C_{15}H_{20}N_5O_{13}P_3S$ (free acid)
MW	603.3 (free acid)
Purity	> 95% HPLC
Quantity	5 μ mol \approx 3mg
Solubility	Sodium salt soluble in water.
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