

p70 S6 kinase (human), (recombinant)

p70 S6K was first identified as the kinase responsible for the phosphorylation of the S6 ribosomal protein. Research has raised the possibility that S6K might, in fact, be the rapamycin-sensitive effector responsible for directly phosphorylating IRS-1 and thereby inhibiting PI 3-kinase activation. Another study suggests that S6K1 activation may mediate some of the effects of a high-fat diet on inhibiting the insulin-PI 3-kinase signaling pathway³. This makes S6K1 a promising new therapeutic target in the fight against type 2 diabetes.

Ordering Information

[Order Online »](#)

BML-SE345-0005	5µg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name p70 S6K

Application Notes Useful for kinetic and functional studies, phosphorylation of target substrates, drug screening.

Formulation Liquid. In 50mM sodium phosphate, pH 7.0, containing 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, and 25% glycerol.

MW 71 kDa

Source Produced in insect cells. Active, full length p70 S6K. Produced in a baculovirus expression system.

Technical Info / Product Notes Replacement for **ADI-PPK-439**

UniProt ID P23443 (S6K1), Q9UBS0 (S6K2)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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