

Lck (human), (recombinant) (GST- tag)

Lck p56 has been shown to phosphorylate proteins that are part of the T cell receptor signaling cascade such as RasGAP, p56.

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 p56lck

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

Formulation Liquid. In 50mM TRIS-HCl, pH 7.5, containing 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

MW 84 kDa

Source Produced in insect cells. Produced in a baculovirus expression system.

Technical Info / Product Notes Replacement for **ADI-PPK-430**

UniProt ID P06239

Last modified: May 29, 2024



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