

# Lyn B (human), (recombinant) (GST- tag)

## Ordering Information

[Order Online »](#)

BML-SE385-0005	5µg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Use/Stability	After initial defrost, aliquot product into individual tubes and refreeze the remaining, unused enzyme quickly by snap-freezing in a dry/ice ethanol bath or liquid nitrogen, if possible.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Tyrosine protein kinase Lyn B
Application Notes	Useful for kinetic and functional studies, phosphorylation of target substrates, drug screening.
Formulation	Liquid. 0.1 mg/ml in 50 mM TRIS-HCl, pH 7.5, containing 150mM sodium chloride, 0.25mM DTT, 0.1mM EGTA, 0.1mM EDTA, 0.1mM PMSF and 25% glycerol.
MW	~85 kDa
Purity	≥90% (SDS-PAGE)
Source	Produced in insect cells. Full length active enzyme. Produced in a baculovirus expression system.
Specific Activity	180 nmol/min/mg. Kinase activity was measured as the molar amount of phosphate incorporated into poly (Glu-Tyr) at 30°C using a concentration of 50µM [ <sup>32</sup> P]ATP in 25mM MOPS, pH 7.2, 12.5mM β-glycerophosphate, 5mM EGTA, 2mM EDTA, 25mM MgCl <sub>2</sub> , 0.25mM DTT (added right before using).
Technical Info / Product Notes	Replacement for <b>ADI-PPK-431</b>
UniProt ID	P07948-2



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-  
be@enzolifesciences.com](mailto:info-be@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)