

PRODUCT INFORMATION



Recombinant Prokaryotic Lectin- α Man (*P. aeruginosa*)

Item No. 36395

Overview and Properties

Synonym:	RPL- α Man
Source:	Recombinant His-tagged <i>P. aeruginosa</i> RPL- α Man expressed in <i>E. coli</i>
Molecular Weight:	13.13 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥95% estimated by SDS-PAGE
Supplied in:	PBS, pH 7.2

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Recombinant prokaryotic lectin- α Man (RPL- α Man) is a *P. aeruginosa* lectin expressed in and purified from *E. coli*. It binds to glycoproteins containing Lewis a (Le^a), Lewis X (Le^X), or terminal α -mannose groups. RPL- α Man activity requires metal ions, therefore, all buffers must be supplemented with 1 mM calcium chloride, magnesium chloride, and manganese chloride. It is recommended to use metal ion-supplemented Tris-buffered saline (TBS), pH 7.6, to dilute RPL- α Man to the desired working concentration (≤ 20 μ g/ml). Following dilution, incubate RPL- α Man for a minimum of 30 minutes prior to use to allow metal complexation.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/29/2022

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM