

# PRODUCT INFORMATION



## Recombinant Prokaryotic Lectin-Man2 (*B. cenocepacia*)

Item No. 35585

### Overview and Properties

<b>Synonym:</b>	RPL-Man2
<b>Source:</b>	Recombinant His-tagged <i>B. cenocepacia</i> RPL-Man2 expressed in <i>E. coli</i>
<b>Molecular Weight:</b>	15.163 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	≥95% estimated by SDS-PAGE
<b>Supplied in:</b>	PBS, pH 7.2

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

### Description

Recombinant prokaryotic lectin-Man2 (RPL-Man2) is a *B. cenocepacia* lectin expressed in and purified from *E. coli*.<sup>1</sup> It is a monomer and selectively binds to glycoproteins containing terminal  $\alpha$ -mannose groups, such as glucose oxidase, over glycoproteins containing  $\alpha(1\rightarrow6)$  fucose or  $\beta(1\rightarrow4)$  galactose groups. RPL-Man2 has been used to detect glycan structures in biolayer interferometry experiments. RPL-Man2 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-Man2 to the desired working concentration ( $\leq 20$   $\mu\text{g/ml}$ ). Following dilution, incubate RPL-Man2 for a minimum of 30 minutes prior to use to allow metal complexation.

### Reference

1. Fernandez-Poza, S., Padros, A., Thompson, R., *et al.* Tailor-made recombinant prokaryotic lectins for characterisation of glycoproteins. *Anal. Chim. Acta* **1155**, 338352 (2021).

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, 05/19/2022

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