

PRODUCT INFORMATION

Anthrax Lethal Factor (recombinant from *Bacillus anthracis*) Modified Sequence, (LF-HMA)

Item No. 23216

Overview and Properties

Contents:	Each vial, when reconstituted with 0.5 ml of water, contains 1 mg of lethal factor from <i>Bacillus anthracis</i> in 5 mM HEPES, 50 mM sodium chloride, pH 7.5.
Storage:	-20°C (as supplied); -20°C or -80°C (after reconstitution); avoid freeze/thaw cycles
Stability:	≥2 years
Purity:	≥90% (estimated by SDS-PAGE)
Protein	
Concentration:	The protein concentration prior to lyophilization was determined by absorbance at 280 nm using an extinction coefficient of 0.88 mg/ml ⁻¹ .

Special Conditions: Handle gently; do not vortex

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

Description

Anthrax lethal factor is one component of a tripartite protein toxin secreted into host systems during anthrax infection caused by *B. anthracis*.¹ The toxin is a mixture of three protein components: Lethal factor, edema factor, and protective antigen. Lethal factor is a zinc-dependent metalloprotease that disrupts MAPK signaling by cleaving MAPK kinases near the N-terminal end. It requires protective antigen, but not edema factor, for its lethality.

Reference

1. Pannifer, A.D., Wong, T.Y., Schwarzenbacher, R., *et al.* Crystal structure of the anthrax lethal factor. *Nature* **414**(6860), 229-233 (2001).

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, 09/24/2019

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