

# PRODUCT INFORMATION



## Extracellular Death Factor

Item No. 15478

<b>CAS Registry No.:</b>	960129-66-2	
<b>Formal Name:</b>	L-asparaginy-L-asparaginy-L-tryptophyl-L-asparaginy-L-asparagine	
<b>Synonym:</b>	EDF	
<b>MF:</b>	$C_{27}H_{36}N_{10}O_{10}$	H—Asn—Asn—Trp—Asn—Asn—OH
<b>FW:</b>	660.6	
<b>Purity:</b>	≥95%	
<b>Supplied as:</b>	A solid	
<b>Storage:</b>	-20°C	
<b>Stability:</b>	≥4 years	

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

### Laboratory Procedures

Extracellular death factor (EDF) is supplied as a solid. A stock solution may be made by dissolving the EDF in the solvent of choice. EDF is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. EDF is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

EDF is a linear pentapeptide that communicating cells produce and release, which upon reaching a sufficient concentration activates the cell death pathway in a subset of cells.<sup>1</sup> It is sensitive to extreme pH, high temperatures, and other stressful conditions.<sup>1</sup> At 2.5 ng/ml, EDF has been shown to facilitate *mazEF*-mediated cell death, significantly reducing population size in *E. coli* cultures.<sup>1</sup>

### Reference

1. Kolodkin-Gal, I., Hazan, R., Gaathon, A., *et al.* A linear pentapeptide is a quorum-sensing factor required for *mazEF*-mediated cell death in *Escherichia coli*. *Science* **318(5850)**, 652-655 (2007).

#### 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, 11/28/2022

#### 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