

# PRODUCT INFORMATION



## Flagellin 22

Item No. 42316

CAS Registry No.: 304642-91-9

Formal Name: L-glutaminy-L-arginyl-L-leucyl-L-seryl-L-threonylglycyl-L-seryl-L-arginyl-L-isoleucyl-L-asparaginy-L-seryl-L-alanyl-L-lysyl-L- $\alpha$ -aspartyl-L- $\alpha$ -aspartyl-L-alanyl-L-alanylglycyl-L-leucyl-L-glutaminy-L-isoleucyl-L-Alanine

H—Gln—Arg—Leu—Ser—Thr—Gly—Ser—Arg—Ile—Asn—  
Ser—Ala—Lys—Asp—Asp—Ala—Ala—Gly—Leu—Gln—  
Ile—Ala—OH

Synonym: Flg22

Peptide Sequence: QRLSTGSRINSAKDDAAGLQIA-OH

MF:  $C_{93}H_{162}N_{32}O_{34}$

FW: 2,272.5

Purity:  $\geq 90\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years

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

### Laboratory Procedures

Flagellin 22 is supplied as a solid. Aqueous solutions of flagellin 22 can be prepared by directly dissolving the solid in aqueous buffers. Flagellin 22 is soluble ( $\geq 10$  mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Flagellin 22 is a peptide pathogen-associated molecular pattern (PAMP) in plants.<sup>1</sup> It inhibits growth, reduces fresh weight, and increases callose levels in *A. thaliana* seedlings when used at a concentration of 10  $\mu\text{M}$ .

### Reference

1. Gómez-Gómez, L., Felix, G., and Boller, T. A single locus determines sensitivity to bacterial flagellin in *Arabidopsis thaliana*. *Plant J.* **18(3)**, 277-284 (1999).

#### 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, 09/04/2025

#### 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