

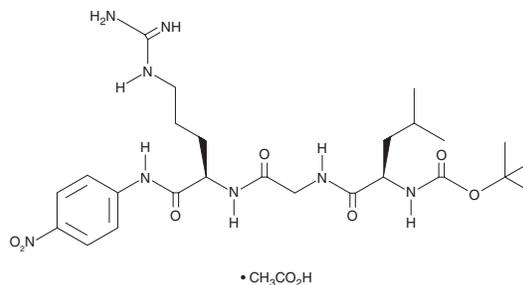
# PRODUCT INFORMATION



## Boc-LGR-pNA (acetate)

Item No. 35176

**CAS Registry No.:** 176786-91-7  
**Formal Name:** N-[(1,1-dimethylethoxy)carbonyl]-L-leucylglycyl-L-(4-nitrophenyl)-L-argininamide, monoacetate  
**Synonyms:** Boc-Leu-Gly-Arg-pNA,  
Boc-Leu-Gly-Arg-p-Nitroanilide,  
Tertbutyloxycarbonyl-Leu-Gly-Arg-p-Nitroanilide  
**MF:** C<sub>25</sub>H<sub>40</sub>N<sub>8</sub>O<sub>7</sub> • C<sub>2</sub>H<sub>4</sub>O<sub>2</sub>  
**FW:** 624.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 202, 222, 312 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Boc-LGR-pNA (acetate) is supplied as a solid. A stock solution may be made by dissolving the Boc-LGR-pNA (acetate) in the solvent of choice, which should be purged with an inert gas. Boc-LGR-pNA (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Boc-LGR-pNA (acetate) in ethanol and DMSO is approximately 15 mg/ml and approximately 10 mg/ml in DMF.

### Description

Boc-LGR-pNA is a chromogenic substrate for the detection of bacterial endotoxins.<sup>1</sup> Upon enzymatic cleavage by endotoxin-induced amidases, p-nitroanilide (pNA) is released which can be quantified by colorimetric detection at 405 nm as a measure of endotoxin levels.

### Reference

1. Harada-Suzuki, T., Morita, T., Iwanaga, S., *et al.* Further studies on the chromogenic substrate assay method for bacterial endotoxins using horseshoe crab (*Tachypleus tridentatus*) hemocyte lysate. *J. Biochem.* **92**(3), 793-800 (1982).

#### 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, 12/12/2022

#### 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