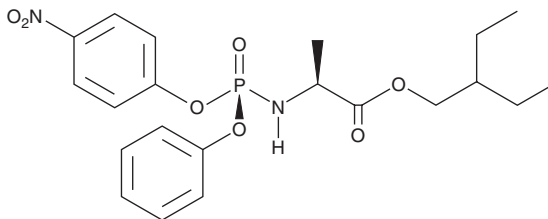


# PRODUCT INFORMATION

## N-[(S)-(4-Nitrophenoxy)phenoxyphosphinyl]-L-alanine 2-ethylbutyl ester Item No. 34748

CAS Registry No.: 1354823-36-1  
MF:  $C_{21}H_{27}N_2O_7P$   
FW: 450.4  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{max}$ : 270 nm  
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

N-[(S)-(4-Nitrophenoxy)phenoxyphosphinyl]-L-alanine 2-ethylbutyl ester is supplied as a solid. A stock solution may be made by dissolving the N-[(S)-(4-nitrophenoxy)phenoxyphosphinyl]-L-alanine 2-ethylbutyl ester in the solvent of choice, which should be purged with an inert gas. N-[(S)-(4-Nitrophenoxy)phenoxyphosphinyl]-L-alanine 2-ethylbutyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-[(S)-(4-nitrophenoxy)phenoxyphosphinyl]-L-alanine 2-ethylbutyl ester in ethanol is approximately 25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

N-[(S)-(4-Nitrophenoxy)phenoxyphosphinyl]-L-alanine 2-ethylbutyl ester is a synthetic intermediate.<sup>1</sup> It has been used in the synthesis of remdesivir (Item No. 30354).

### Reference

- Warren, T.K., Jordan, R., Lo, M.K., *et al.* Therapeutic efficacy of the small molecule GS-5734 against Ebola virus in rhesus monkeys. *Nature* **531**(7594), 381-385 (2016).

#### 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, 11/10/2022

#### 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