

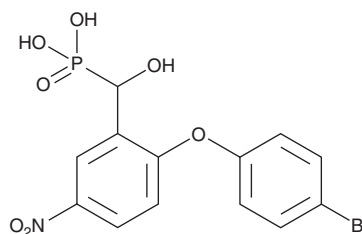
# PRODUCT INFORMATION



**RWJ 60475**

Item No. 45483

**CAS Registry No.:** 204130-08-5  
**Formal Name:** P-[[2-(4-bromophenoxy)-5-nitrophenyl]hydroxymethyl]-phosphonic acid  
**MF:** C<sub>13</sub>H<sub>11</sub>BrNO<sub>7</sub>P  
**FW:** 404.1  
**Purity:** ≥98%  
**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

RWJ 60475 is supplied as a solid. A stock solution may be made by dissolving the RWJ 60475 in the solvent of choice, which should be purged with an inert gas. RWJ 60475 is sparingly soluble (1-10 mg/ml) in methanol and DMSO.

## Description

RWJ 60475 is an inhibitor of CD45 (IC<sub>50</sub> = 2 μM).<sup>1</sup>

## Reference

1. Beers, S.A., Malloy, E.A., Wu, W., *et al.* Nitroarylhydroxymethylphosphonic acids as inhibitors of CD45. *Bioorg. Med. Chem.* **5(12)**, 2203-2211 (1997).

### 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, 04/09/2026

## 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