

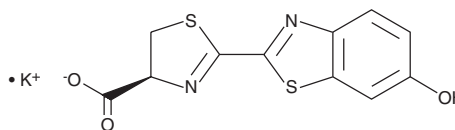
PRODUCT INFORMATION



D-Luciferin (potassium salt)

Item No. 14681

CAS Registry No.: 115144-35-9
Formal Name: 4,5-dihydro-2-(6-hydroxy-2-benzothiazolyl)-4S-thiazolecarboxylic acid, monopotassium salt
MF: C₁₁H₇N₂O₃S₂ • K
FW: 318.4
Purity: ≥95%
UV/Vis.: λ_{max}: 268, 327 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

D-Luciferin (potassium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the D-luciferin (potassium salt) in the solvent of choice. D-Luciferin (potassium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of D-luciferin (potassium salt) in these solvents is approximately 0.25, 10, and 16.7 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of D-luciferin (potassium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of D-luciferin (potassium salt) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

D-Luciferin is a chemiluminescent substrate of firefly luciferase.¹ It produces light upon oxidative decarboxylation by luciferase in the presence of ATP. D-Luciferin can be employed to assay the expression of the luciferase gene linked to a promoter of interest. Alternatively, D-luciferin and luciferase can be used to assess ATP availability in cellular or biochemical assays.

Reference

1. Lember, N. Firefly luciferase can use L-luciferin to produce light. *Biochem. J.* **317(1)**, 273-277 (1996).

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, 09/25/2018

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