

PRODUCT INFORMATION



Luciferase (firefly, recombinant)

Item No. 20398

CAS Registry No.: 61970-00-1

Purity: ≥95% estimated by SDS-PAGE

Molecular Weight: 60 kDa

Supplied as: A 1 mg/ml solution in 50 mM Tris buffer with 30% glycerol, pH 7.5

Storage: -20°C

Stability: ≥2 years

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

Description

Luciferase (firefly, recombinant) is a light-generating enzyme that catalyzes an oxidative reaction, using luciferin, ATP, and molecular oxygen to produce oxyluciferin, which yields visible light.^{1,2} The bioluminescence of this ATP-dependent luciferase is commonly used in reporter systems that are amenable to high-throughput screening.³

References

1. Baldwin, T. O. Firefly luciferase: The structure is known, but the mystery remains. *Structure* **4(3)**, 223-228 (1996).
2. Xu, T., Close, D., Handagama, W., *et al.* The expanding toolbox of *in vivo* bioluminescent imaging. *Front. Oncol.* **6**, 150 (2016).
3. Auld, D. S. and Ingles, J. Interferences with luciferase reporter enzymes. *Assay Guidance Manual* (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, 11/02/2022

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