

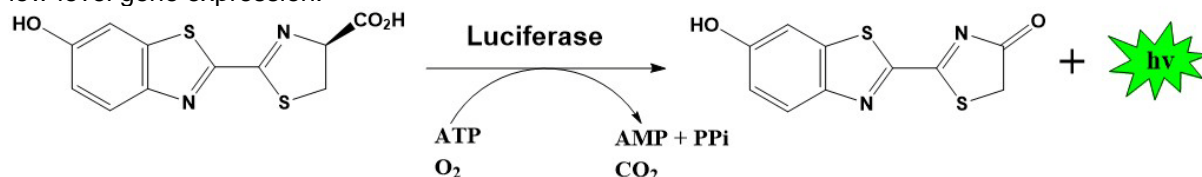
## D-Luciferin

Catalog Number	Product Name	Packaging Size
C292A	D-Luciferin, free acid	10 mg
C292B	D-Luciferin, free acid	25 mg
C292C	D-Luciferin, free acid	100 mg
C293A	D-Luciferin, potassium salt	10 mg
C293B	D-Luciferin, potassium salt	25 mg
C293C	D-Luciferin, potassium salt	100 mg
C294A	D-Luciferin, sodium salt	10 mg
C294B	D-Luciferin, sodium salt	25 mg
C294C	D-Luciferin, sodium salt	100 mg

**Storage upon receipt:** -20°C, protected from light

### Introduction

**Luciferins** are a class of ATP-dependent substrates that are oxidized in the presence of the enzyme luciferase to produce oxyluciferin and energy in the form of light. Luciferin undergoes an enzyme-catalysed oxidation and the resulting unstable reaction intermediate emits light upon decaying to its ground state. This system is employed as a very useful reporter in plants, bacteria, and mammalian cells. Because chemiluminescent techniques are virtually background-free, this reporter gene system is ideal for detecting low-level gene expression.



### Specifications

Product Name:	D-Luciferin, free acid	D-Luciferin, potassium salt	D-Luciferin, sodium salt
Molecular Formula:	C <sub>11</sub> H <sub>8</sub> N <sub>2</sub> O <sub>3</sub> S <sub>2</sub>	C <sub>11</sub> H <sub>7</sub> KN <sub>2</sub> O <sub>3</sub> S <sub>2</sub>	C <sub>11</sub> H <sub>7</sub> N <sub>2</sub> NaO <sub>3</sub> S <sub>2</sub>
Molecular Weight:	280.32	318.40	302.30
CAS Number:	2951-17-5	15144-35-9	103404-75-7
Storage Conditions:	-20 °C, protected from light	-20 °C, protected from light	-20 °C, protected from light
Shipping Condition:	Room Temperature	Room Temperature	Room Temperature
Structure:			

## References:

1. Bacterial and Firefly Luciferase Genes in Transgenic Plants, Advantages and Disadvantages of a Reporter Gene.  
Koncz C, et al.  
Dev Genet (1990) 11:224-224
2. Investigation of the Interaction between Firefly Luciferase and Oxyluciferin or Its Analogues by Steady State and Subnanosecond Time-Resolved Fluorescence. Investigation of the Interaction between Firefly Luciferase and Oxyluciferin or Its Analogues by Steady State and Subnanosecond Time-Resolved Fluorescence.  
Gandelman OA, et al.  
J Photochem Photobiol B (1994) 22:203-203
3. ATP Determination with Firefly Luciferase. ATP Determination with Firefly Luciferase.  
Leach FR  
J Appl Biochem (1981) 3:473-473