

**D-Luciferin phosphate \*CAS 145613-12-3\***

Catalog number: 12512

Unit size: 1 mg

Component	Storage	Amount
D-Luciferin phosphate *CAS 145613-12-3*	Freeze (<-15 °C), Minimize light exposure	1 mg

**OVERVIEW**

D-Luciferin phosphate is a highly sensitive bioluminescent substrate for detecting alkaline phosphatases and protein phosphatases by luminometric detection. It is used in combination with luciferase/ATP system. D-Luciferin phosphate can be converted to D-Luciferin by phosphatases (such as alkaline phosphatases, acid phosphatases and protein phosphatases) that gives luminescence in the presence of luciferase and its cofactors. Our Amplite luciferase assay kit can be used for detecting D-luciferin release from D-Luciferin phosphate by phosphatases.

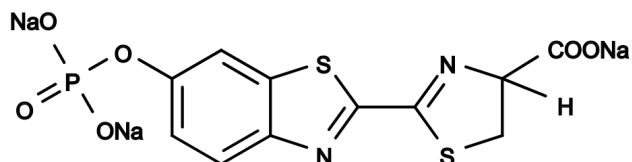
**AT A GLANCE**

**Important** It is important to store at <-15 °C and should be stored in cool, dark place.

It can be used within 12 months from the date of receipt.

**PREPARATION OF WORKING SOLUTION**

Working solution should be prepared in water or DMSO.

**EXAMPLE DATA ANALYSIS AND FIGURES**


**Figure 1.** Chemical structure for D-Luciferin phosphate \*CAS 145613-12-3\*

**DISCLAIMER**

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.