

## **D-Luciferin acetate**

Catalog number: 12516

Unit size: 5 mg

## **Product Details**

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

## **Chemical Properties**

Appearance White solid

Molecular Weight 322.36

Soluble In DMSO

**Chemical Structure** 

## **Applications**

D-Luciferin acetate contains an aceate moiety attached at the 6-O-position, thus preventing it be recognized by the firefly luciferase enzyme. However, the acatate moiety is removed effectively by esterase activity, and the resulted D-luciferin is well recognized by luciferase. D-Luciferin acetate is a sensitive substrate for the chemiluminescent measurement of esterase activity in homogeneous assays, or in cell lysate samples in combination with luciferase and its cofactors.