

Product Information Sheet

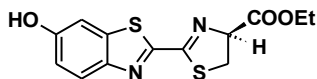
Ordering Information

Product Number: 12515
Product Name: D-Luciferin ethyl ester *Single isomer*
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 308.38

Chemical Structure:



Soluble in: DMSO

Application Notes

D-Luciferin ethyl ester contains an ethyl moiety attached at the carboxy position, thus preventing it be recognized by the firefly luciferase enzyme. However, the ester is hydrolyzed effectively by lipase activity, and the resulted D-luciferin is well recognized by luciferase. D-Luciferin ethyl ester is a sensitive substrate for the chemiluminescent measurement of lipase activity in homogeneous assays, or in cell lysate samples in combination with luciferase and its cofactors.