

Recombinant Obelin

Catalog Number L001-50UG

Catalog Number L001-100UG

Catalog Number L001-250UG



ARBOR
ASSAYS

FEATURES

- Photoprotein from the jellyfish, *Obelia longissima*
- Recombinant protein expressed in *E. coli*
- High specific activity

INTRODUCTION

Recombinant bioluminescent photoprotein from *Obelia longissima*. The photoprotein is approximately 21,000 molecular weight and contains the luminescent substrate, native coelenterazine, and oxygen bound to the protein. Upon addition of excess calcium ions that fill the three calcium binding sites, the oxygen reacts with the substrate to yield a rapid flash of blue light centered at 485 nm. This emission reaches a peak within 100 msec. and decays in less than 1 second.

FORM:	Lyophilized protein from Lyophilization Buffer.
BUFFER COMPOSITION:	10 mM Tris, 1 mM EDTA, 0.1% Tween 20, 10% mannitol, pH 7.5.
PURITY:	Purified to homogeneity.
ACTIVITY:	$> 0.5 \times 10^{12}$ RLU/mg*
STORAGE:	4°C for short term, up to 24 months at -20°C.
USES:	Calcium ion determination.
COUNTRY OF ORIGIN:	USA

* As measured for 1 second on a Berthold Orion II plate luminoter using 150 μ L of Arbor Assays Obelin Trigger Solution, Catalog Number X006- 15ML or X006-100ML.

FOR RESEARCH USE ONLY