

Highly Stable ATP Assay Kit ab287863

2 Images

Overview

Product name	Highly Stable ATP Assay Kit
Detection method	Luminescent
Sample type	Tissue
Assay type	Enzyme activity
Sensitivity	> 1 nM
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mammals
Product overview	<p>In this kit, Highly Stable ATP Assay Kit ab287863 (previously known as StayBrite Higly stable ATP Bioluminescence assay kit, K791), we have developed a highly stable Luciferase, which is a genetically modified variant derived from the Luciferase of Diaphanes pectinealis (Chinese Firefly) endemic to Yunnan province, China.</p> <p>We named the highly stable luciferase rLucHS. Compared to the normal phenotype of Photinus pyralis, the rLucHS provides enhanced stability, excellent sensitivity, and a broader and more physiologically relevant effective pH range. At all pH's below ~8.2, rLucHS is significantly more active than Photinus luciferase and is stable for weeks at room temperature and >60 minutes at 37°C. The activity profile shows a linear detection range for ATP values as low as 10 fmol/assay (1 nM). The specific activity of rLucHS is ~ 5x10⁴ RLU/mg protein. The assay can be fully automated for high throughput (1 second/sample) and is extremely sensitive. The high sensitivity of this assay is ideal for detecting ATP production or consumption in various enzymatic reactions.</p>
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K791 StayBrite™ Highly Stable ATP Bioluminescence Assay Kit. K791-100 is the same size as the 100 test size of ab287863.</p>
Platform	Microplate (12 x 8 well strips)

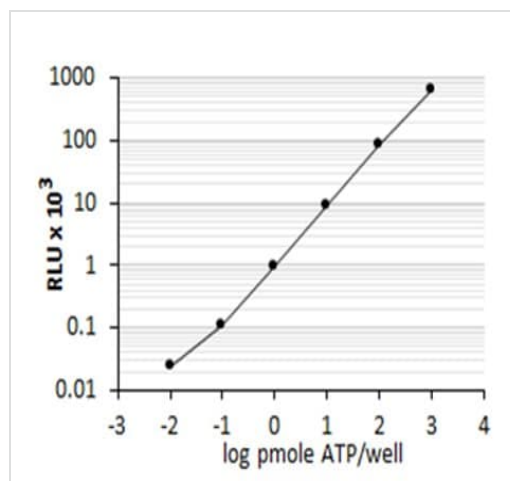
Properties

Storage instructions	Store at -20°C. Please refer to protocols.
----------------------	--

Components	1000 tests
10X StayBite Reaction buffer	1 x 20ml

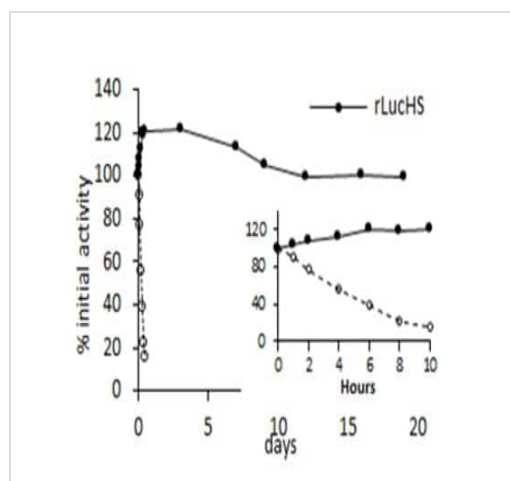
Components	1000 tests
ATP Standard (MW 551) (Lyophilized)	1 vial
StayBite Enzyme Mix (Lyophilized)	1 vial
StayBite reconstitution buffer	1 x 11ml

Images



Standard curve analysis of ab287863. For reference use only.

Highly Stable ATP Assay Kit - Standard Curve



rLucHS Stability at Room Temperature.

Highly Stable ATP Assay Kit

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet

- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors