

Renin Assay Kit (Fluorometric) ab138875

4 References 1 图像

概述									
产品名称	Renin Assay试剂盒(Fluorometric)								
检测方法	Fluorescent								
样品类型	Cell Lysate								
检测类型	Quantitative								
灵敏度	10 ng/ml								
种属反应性	与反应: Mammals, Other species								
产品概述	Renin Assay Kit (Fluorometric) (ab138875) provides a convenient assay for high throughput screening of renin inhibitors and for continuous assay of renin activity using a TF3/TQ3 FRET peptide. In the FRET peptide the fluorescence of TF3 is quenched by TQ3. Upon cleavage into two separate fragments by renin, the fluorescence of TF3 is recovered, the fluorescent signal can be easily monitored by a fluorescence microplate reader at Ex/Em = 540/590 nm. This assay is about fifty fold more sensitive than an EDANS/DABCYL-based assay.								
平台	Microplate reader								
性能									
存放说明	Store at -20°C. Please refer to protocols.								
<table><tr><th>组件</th><th>100 tests</th></tr><tr><td>Assay Buffer</td><td>1 x 10ml</td></tr><tr><td>Renin Red Substrate (100x)</td><td>1 x 50µl</td></tr><tr><td>Renin Standard</td><td>1 x 25µl</td></tr></table>		组件	100 tests	Assay Buffer	1 x 10ml	Renin Red Substrate (100x)	1 x 50µl	Renin Standard	1 x 25µl
组件	100 tests								
Assay Buffer	1 x 10ml								
Renin Red Substrate (100x)	1 x 50µl								
Renin Standard	1 x 25µl								
功能	Renin is a highly specific endopeptidase, whose only known function is to generate angiotensin I from angiotensinogen in the plasma, initiating a cascade of reactions that produce an elevation of blood pressure and increased sodium retention by the kidney.								
疾病相关	Defects in REN are a cause of renal tubular dysgenesis (RTD) [MIM:267430]. RTD is an autosomal recessive severe disorder of renal tubular development characterized by persistent fetal anuria and perinatal death, probably due to pulmonary hypoplasia from early-onset oligohydramnios (the Potter phenotype).								

Defects in REN are the cause of familial juvenile hyperuricemic nephropathy type 2 (HNFJ2) [MIM:613092]. It is a renal disease characterized by juvenile onset of hyperuricemia, slowly progressive renal failure and anemia.

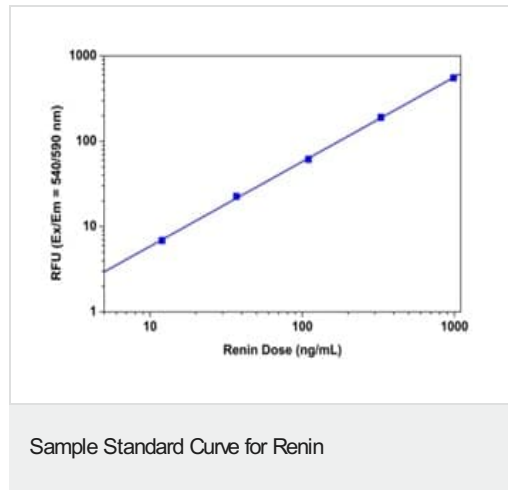
序列相似性

Belongs to the peptidase A1 family.

细胞定位

Secreted. Membrane. Associated to membranes via binding to ATP6AP2.

图片



Renin dose response was measured with ab138875 in a 96-well black solid plate using a fluorescence microplate reader. As low as 10 ng/mL Renin was detected with 60 minutes incubation in 37°C

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.cn/abpromise> or contact our technical team.

Terms and conditions

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