

# ultra-sensitive...

& many other good points!

homogeneous | long wavelengths

as fast as 30 min. | cost effective



## SensoLyte<sup>®</sup> 520 Human Renin Assay Kit

*40x more sensitive*

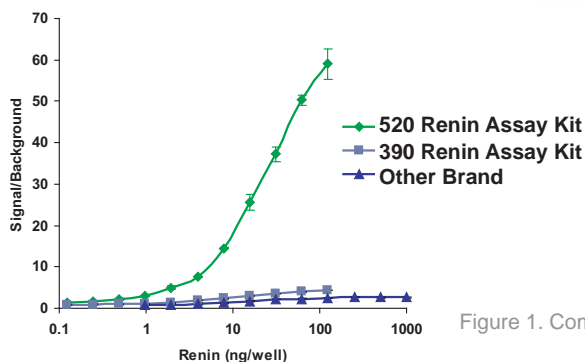


Figure 1. Comparison of 3 renin assay kits.

Sensitivity is a good thing when it comes to high performance assay kits. Offering 40 times greater sensitivity than the next leading brand, AnaSpec's SensoLyte<sup>®</sup> 520 Human Renin Assay Kit is the most sensitive renin assay on the market. Combining high performance with speed and convenience, AnaSpec's series of SensoLyte<sup>®</sup> Renin Assay kits are homogeneous, continuous assays that are optimized for HTS. With a focus on innovation that matches the caliber of our customers, AnaSpec is perfecting the science of trust.



Certified to ISO 9001:2008

# Renin Assay Kits:

Table 1. Comparison data from SensoLyte® 520 & 390 and Company B's renin assay kits.

Characteristics	SensoLyte® 520 Renin Assay Kit, Cat # 72040	SensoLyte® 390 Renin Assay Kit, Cat # 72039	Company B
FRET Pair	5-FAM/QXL™ 520	Dnp/Amp	EDANS/DABCYL
Ex/Em (nm)	490/520	330/390	335/495
Min. Detectable Renin Conc. (ng/ml)	0.8	8	32
Interference from Autofluorescence	Minimal	Moderate	Moderate
Sensitivity	Ultra-sensitive	Good	Minimal

## Related Products:

SensoLyte® 390 ACE2 Activity Assay Kit (Cat# 72086)

Human Renin, Recombinant Protein (Cat # 72041)

# SensoLyte® 520 Rat and Mouse Renin Assay Kits

In addition to the SensoLyte® Human Renin Assay Kits, AnaSpec also provides SensoLyte® 520 Renin Assays specific for rat (Cat# 72140) and mouse (Cat# 72161).

Table 2. SensoLyte® 520 Rat and Mouse renin assay kits.

Characteristics	SensoLyte® 520 Rat Renin Assay Kit Cat # 72140	SensoLyte® 520 Mouse Renin Assay Kit Cat # 72161
FRET Pair	5-FAM/QXL™ 520	5-FAM/QXL™ 520
Ex/Em (nm)	490/520	490/520

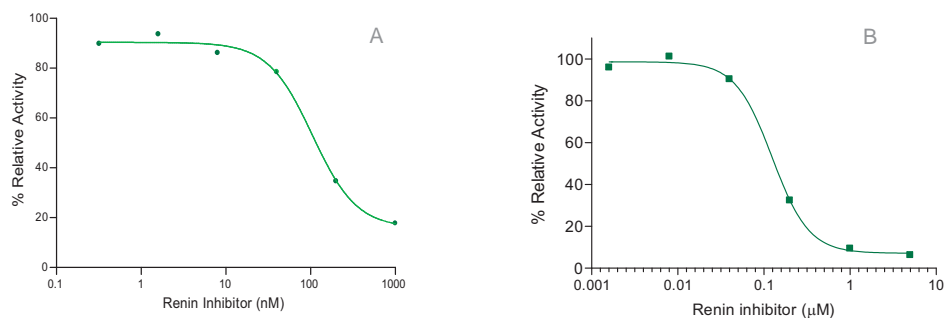


Figure 2. Inhibition results as measured by SensoLyte® 520 Rat (Panel A) and Mouse (Panel B) Renin Assay Kits.

## Related Products:

Rat Renin, Recombinant Protein (Cat # 72141)

Mouse Renin, Recombinant Protein (Cat # 72174)

Rat Renin Inhibitor Peptide, WFML Peptide (Cat # 60463-1)

Angiotensins & Related Peptides

Anti-Renin's (Cat# 54371-54373)