

---

# FGR

Product code: 08-166

Full-length human FGR [1-529(end) amino acids of accession number NP\_005239.1] was expressed as N-terminal GST-fusion protein (86 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP ( $\mu$ M) Km app / Bin	: 34 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.3
IC50 at 1 mM ATP (nM)	: 16

