

---

## AurC(AURKC)

Product code: 05-103

Full-length human AurC [1-275(end) amino acids of accession number NP\_003151.2] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Kemptide
ATP ( $\mu$ M) Km app / Bin	: 24 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.1
IC50 at 1 mM ATP (nM)	: 18

