

---

---

## ARG(ABL2)

Product code: 08-102

Full-length human ARG [2-1161(end) amino acids of accession number NP\_001161708.1] was expressed as N-terminal GST-fusion protein (153 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: ABLtide
ATP ( $\mu$ M) Km app / Bin	: 24 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 27
IC50 at 1 mM ATP (nM)	: 400

