

---

---

# MOS

Product code: 05-118

Full-length, human MOS [1-346(end) amino acids of accession number NP\_005363.1] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system.

Assay platform	:	MSA
Substrate	:	MAP2K1 Cascade Assay *
ATP ( $\mu$ M)	:	1000
Metal	:	Mg
Reference compound	:	Staurosporine
IC50 at 1 mM ATP (nM)	:	32

\* MAP2K1/Erk2/Modified Erktide

---

---