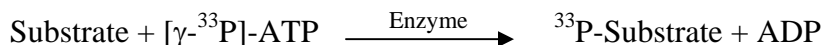


## BRAF

(B-Raf proto-oncogene serine/threonine-protein kinase, B-Raf proto-oncogene serine/threonine-protein, kinase, BRAF1, RAFB1)

- **CAT#:** BRAF
- **Enzyme:** Human BRAF
- **Substrate:** MEK1 (K97R), 1  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
GW 5074	6.2

- **Scientific Information:** P15056 (<http://www.expasy.org/uniprot/P15056>)

### Reference Compound IC50 for BRAF

