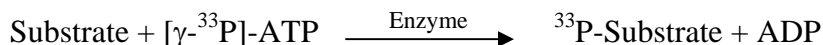


## BRAF (V599E)

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

- **CAT#:** BRAF (V599E)
- **Enzyme:** Human BRAF (V599E)
- **Substrate:** MEK1 (K97R), 1  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds | IC50 (nM) |
|-----------|-----------|
| GW 5074   | 64        |

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

### Reference compound IC50 for BRAF (V599E)

