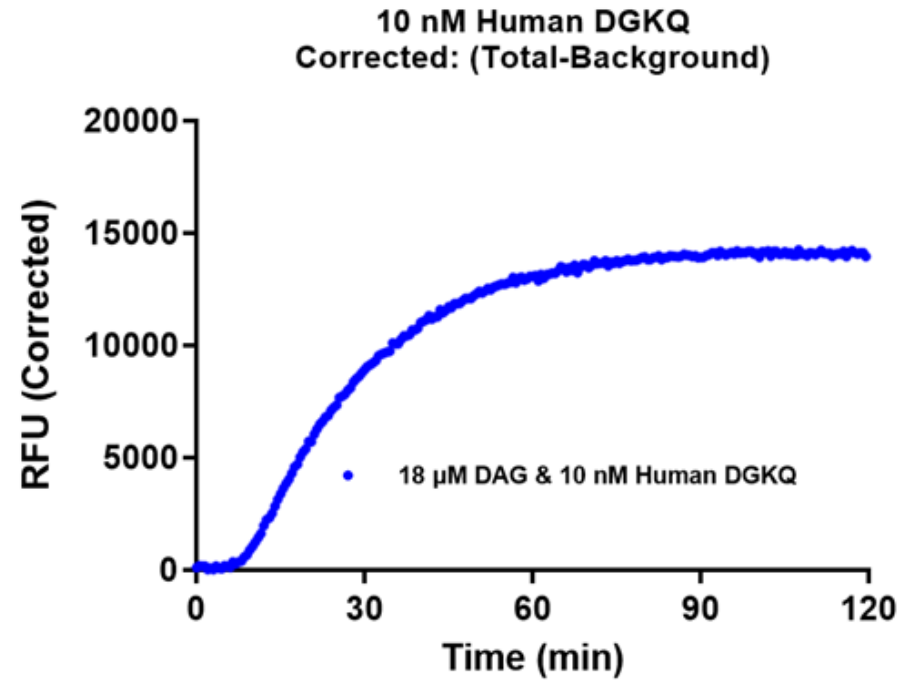
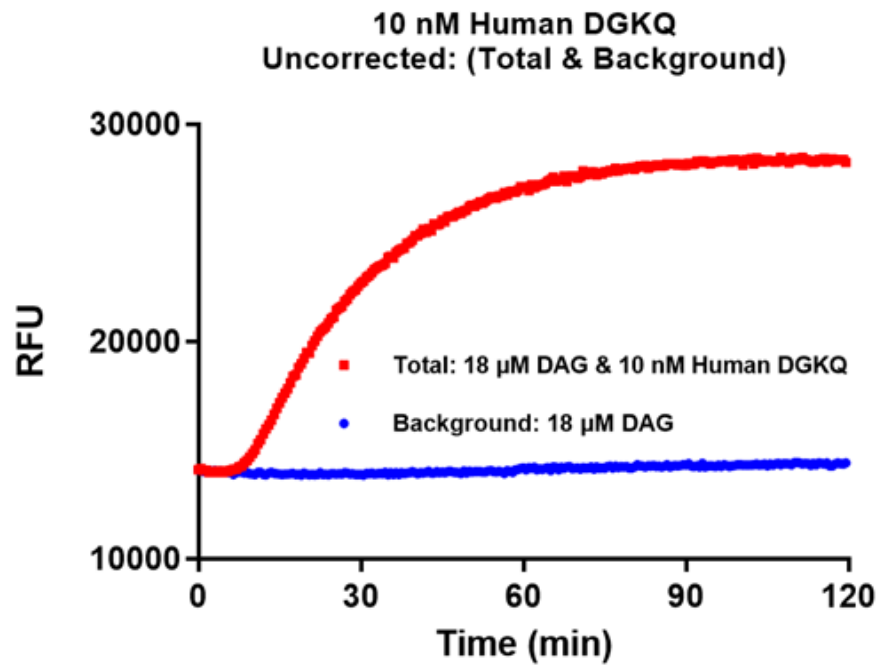


Kinase      BTN-DGKθ(DGKQ)  
Cat.#      12-409-20N  
Substrate   PhosphoSens<sup>®</sup> DAG Substrate (Assay Quant Technologies Inc.)



DAG Reaction rate:  $901 \pm 9$  RFU/pmole/min