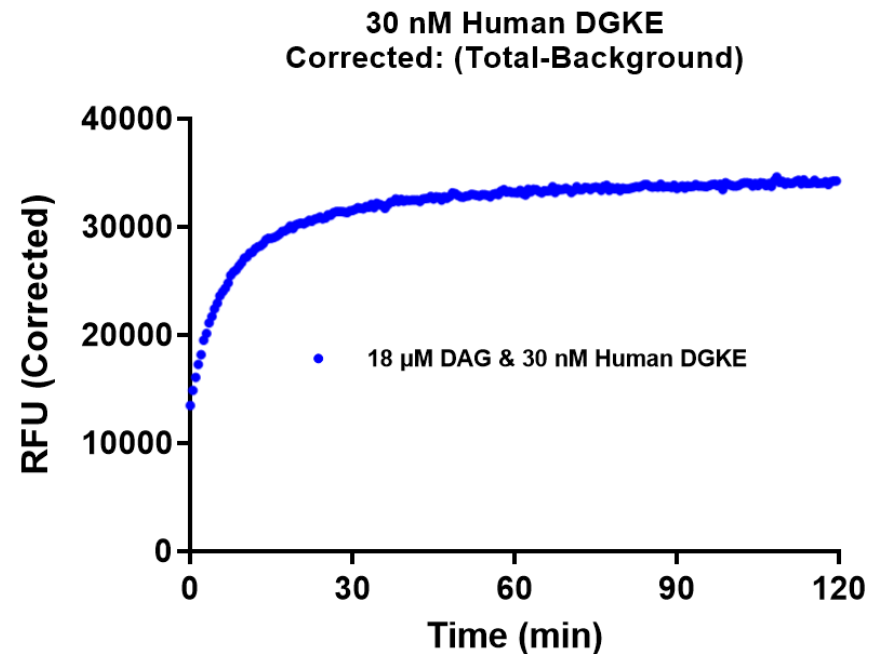
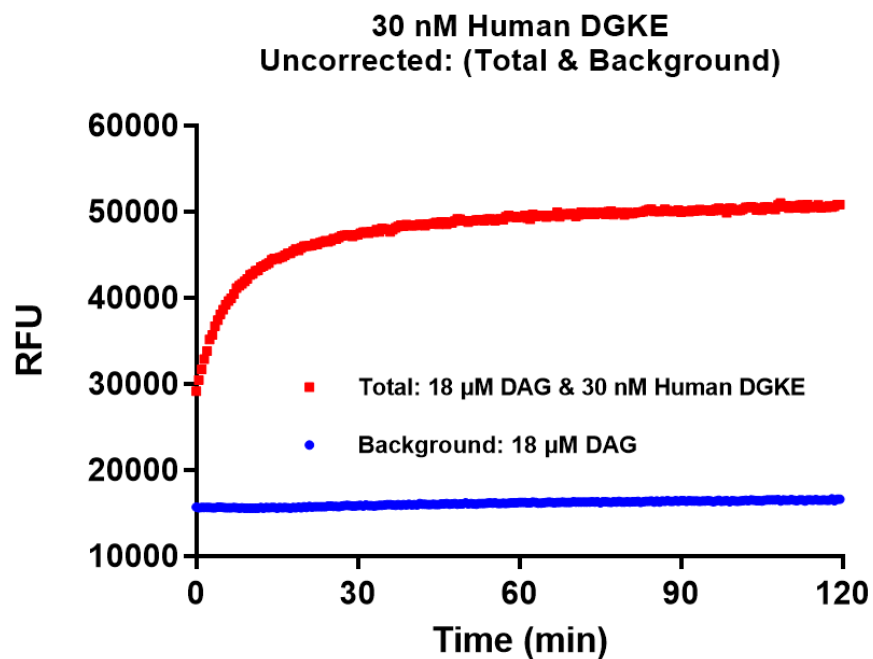


Kinase BTN-DGK ϵ (DGKE)
Cat.# 12-415-20N
Substrate PhosphoSens[®] DAG Substrate (Assay Quant Technologies Inc.)



DAG Reaction rate: $1,261 \pm 56$ RFU/pmole/min