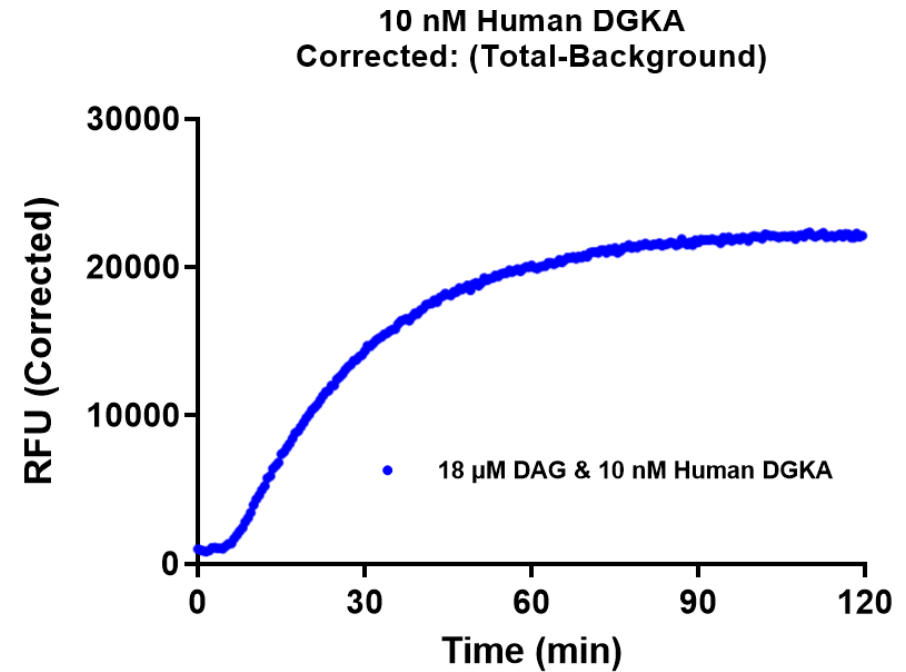
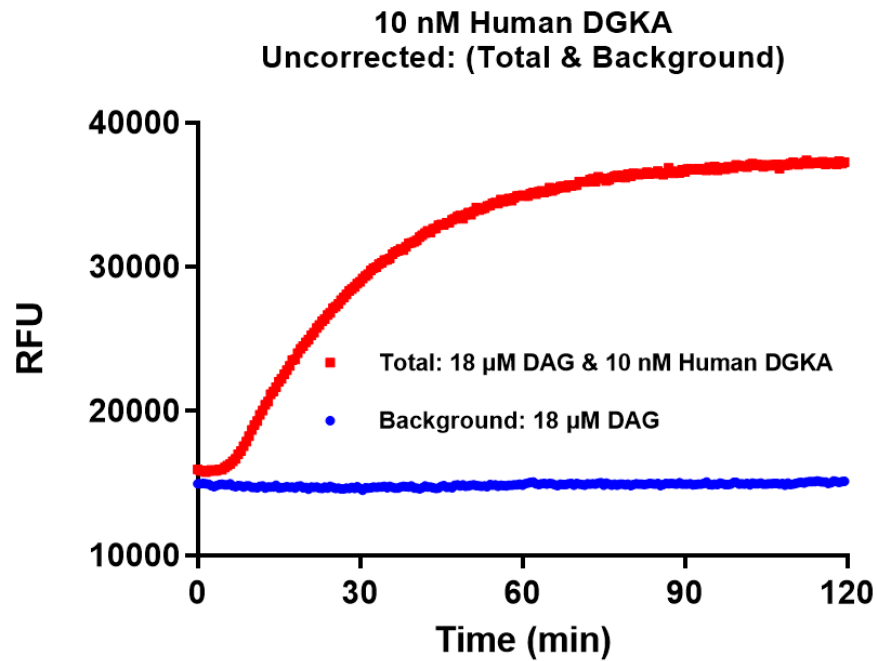


Kinase BTN-DGK α
Cat.# 12-401-20N
Substrate PhosphoSens[®] DAG Substrate (Assay Quant Technologies Inc.)



DAG Reaction rate: $1,230 \pm 17$ RFU/ μ mole/min