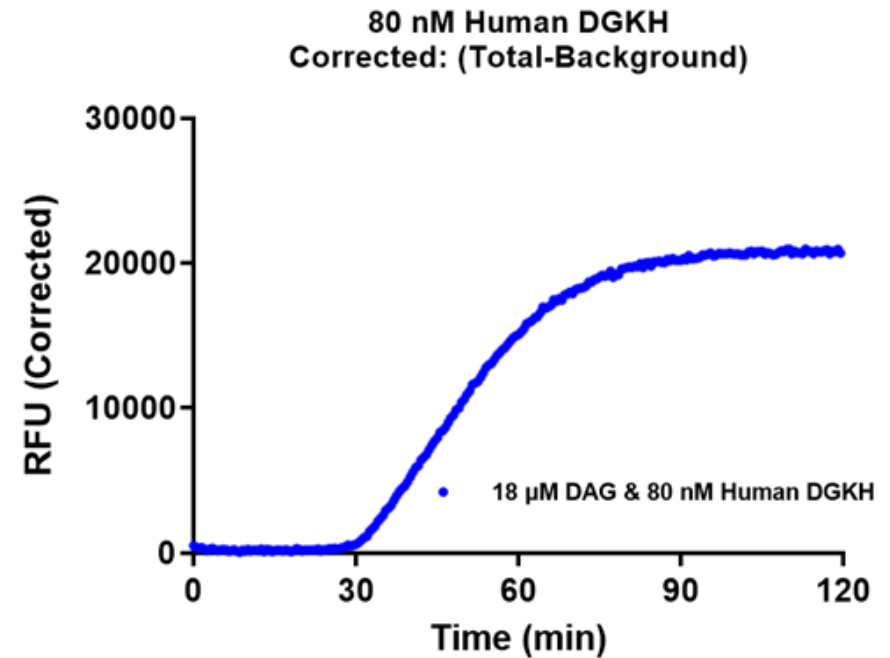
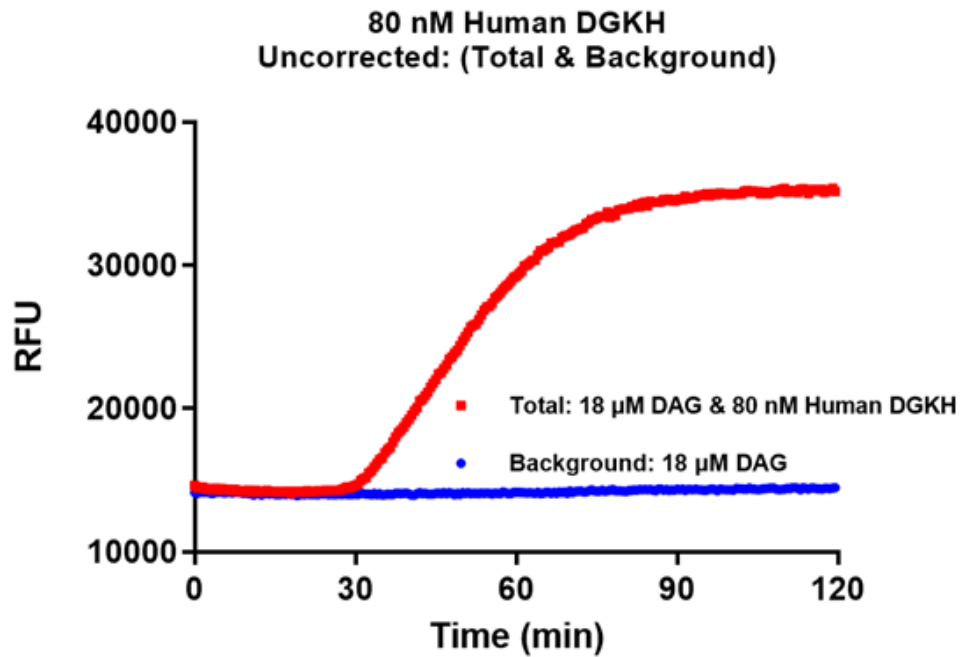


Kinase BTN-DGK η (DGKH)
Cat.# 12-406-20N
Substrate PhosphoSens[®] DAG Substrate (Assay Quant Technologies Inc.)



DAG Reaction rate: 132.9 ± 0.5 RFU/pmole/min