



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Human  $\alpha$ -Synuclein Calibrator  
**Catalog Number:** C01WK-2  
**Lot Number:** A0080167  
**Expiration Date:** 30 SEP 2020  
**Storage Temperature:** -70°C

### Concentration of Materials Tested:

	Assay
	hu $\alpha$ -Synuclein
Calibrator (pg/mL)	208 000

**Concentration:** 20X

**Statement:** This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	24 MAY 2018