



## DECLARATION OF CONFORMITY

(According to ISO/IEC GUIDE 22 and EN 45014)

Manufacturer's Name: *ProteinSimple (a Bio-Techne brand)*

Manufacturer's Address: **3001 Orchard Parkway  
San Jose, CA 95134  
U.S.A.**

### ***DECLARES, THAT THE PRODUCT***

Product Name: *Protein analysis equipment*

Model Number: *Ella*

### ***CONFORMS TO THE FOLLOWING EUROPEAN DIRECTIVES***

***Low Voltage Directive: 2006 / 95 / EC  
EMC Directive: 2004 / 108 / EC  
RoHS Directive: 2011 / 65 / EU***

Supplementary Information:

***Safety: IEC/EN 61010-1:2010; CAN/CSA-C22.2 No. 61010-1:2012***

***EMC: IEC/EN 61326-1:2006***

***Laser Safety: IEC/EN 60825-1:2014***

***Year CE Mark Affixed: 2015***

I, the undersigned, hereby declare that the equipment specified above conforms to the above Directives and Standards.

Place: Wallingford, CT

Signature: \_\_\_\_\_

Date: 08-Sep-2017

Full Name: Henry Couture

Position: Director of Quality