

SelectScience®
The Fastest Way to Expert Opinion



Bronze Seal of Quality

Is hereby granted to

Tecan

for consistently receiving multiple positive reviews on
SelectScience® for the

Tecan Freedom EVO®

A handwritten signature in black ink that reads "K Parker".

Kerry Parker
Editor-in-Chief, SelectScience®



Date awarded: May 30, 2017