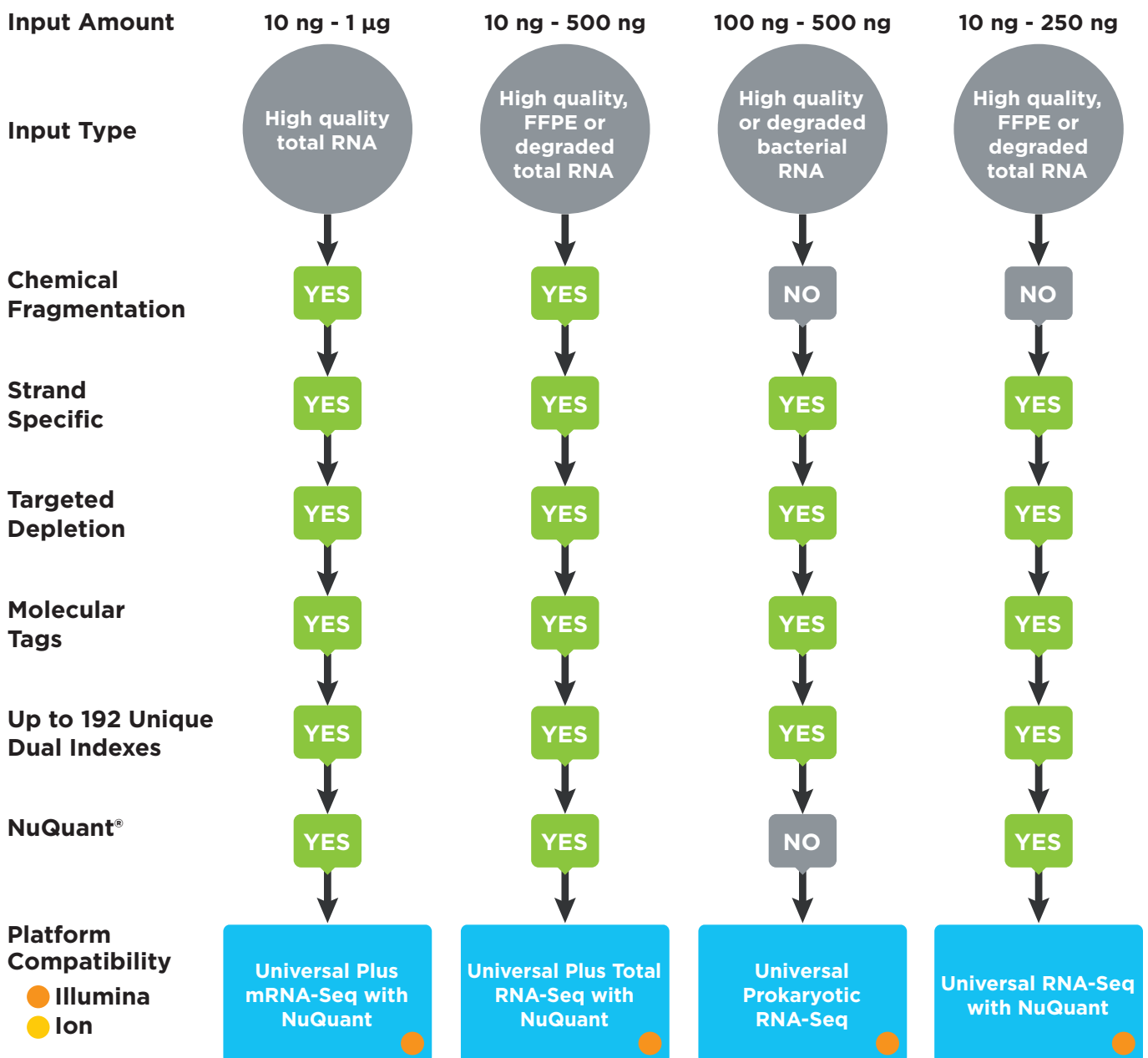


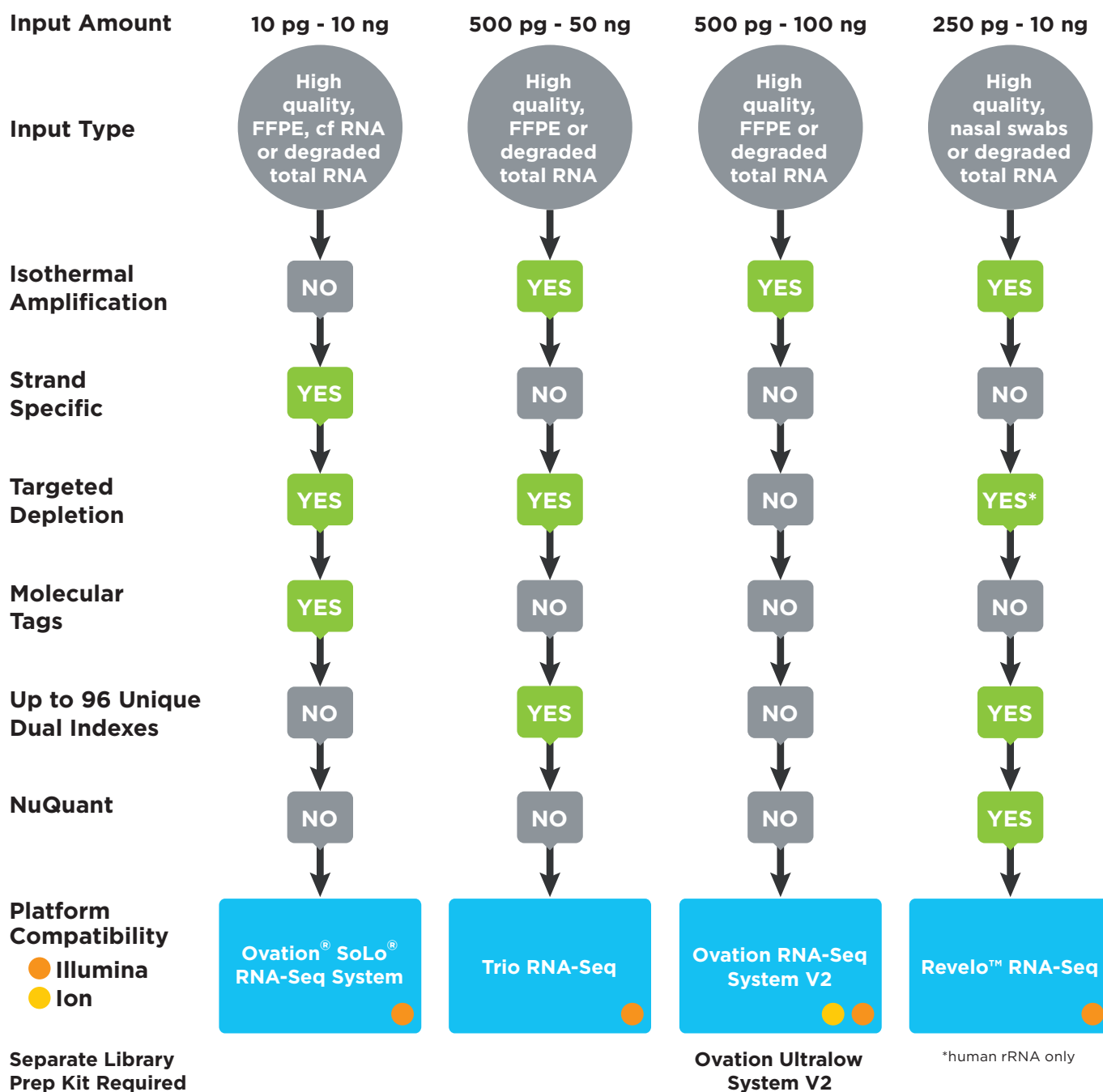
RNA-Seq Solutions.

PRODUCT SELECTION GUIDE

Standard input library preparation solutions for Illumina sequencing



Ultra-low input library preparation solutions for Illumina sequencing



For Research Use Only. Not for use in diagnostic procedures.

www.tecan.com

Australia +61 3 9647 4100 Austria +43 62 46 89 330 Belgium +32 15 42 13 19 China +86 21 220 63 206 France +33 4 72 76 04 80 Germany +49 79 51 94 170 Italy +39 02 92 44 790 Japan +81 44 556 73 11 Netherlands +31 18 34 48 17 4 Nordic +46 8 750 39 40 Singapore +65 644 41 886 Spain +34 93 595 25 31 Switzerland +41 44 922 89 22 UK +44 118 9300 300 USA +1 919 361 5200 Other countries +41 44 922 81 11

Tecan, NuQuant, SoLo, and Ovation are registered trademarks and trademarks of Tecan Group Ltd., Männedorf, Switzerland or of Tecan Genomics, Inc., Redwood City, USA.

© 2020 Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit www.tecan.com

