

IDP-4 Dry Scroll Pump Quick Reference Guide



UNBOXING AND INSTALLATION



Visit the product web page:
[IDP-4 Dry Scroll Pump](#)

This information is subject to change without notice.