

Comparative Specifications

Diagenode sa

CHU, Tour GIGA B34, 3^e étage

Avenue de l'Hôpital, 1

4000 Liège - Belgium

Bioruptor™ UCD-200



Bioruptor™ XL



Adaptors for all types of tubes (0.5, 1.5, 10, 15 and 50ml)	Adaptors for all types of tubes (0.5, 1.5, 10, 15 and 50ml)
Up to 12 tubes of 0.5ml at the same time	Up to 48 tubes of 0.5ml at the same time
Up to 6 tubes of 1.5ml at the same time	Up to 24 tubes of 1.5ml at the same time
Up to 6 tubes of 10-15ml at the same time	Up to 24 tubes of 10-15ml at the same time
Up to 3 tubes of 50ml at the same time	Up to 12 tubes of 50ml at the same time
1 gear plate	2 gear plates
Max sample volume: <ul style="list-style-type: none"> · for 0.5ml tube: 100µl · for 1.5ml tube: 300µl · for 10ml tube: 2ml · for 15ml tube: 2ml · for 50ml tube: 20ml 	Max sample volume: <ul style="list-style-type: none"> · for 0.5ml tube: 100µl · for 1.5ml tube: 300µl · for 10ml tube: 2ml · for 15ml tube: 2ml · for 50ml tube: 20ml
Analogic timer	Digital timer
Plastic noise reduction box	Metal noise reduction box
3 intensity levels (L:160w, M:200w, H:320w)	4 intensity levels (L:130w, M1:200w, M2: 250w, H:300w)
Water cooling by ice or circulation system	Water cooling by ice or circulation system

For more information, please visit our Website www.diagenode.com or :

For Europe and Rest of the World, call +32 4 364 20 50 or send an email to info@diagenode.com

For North America, call +1 973 726 3290 or send an email to infousa@diagenode.com