



BD

BD Mouse CD Marker Chart

bdbiosciences.com/cdmarkers

