

Manual document for all ACD products.

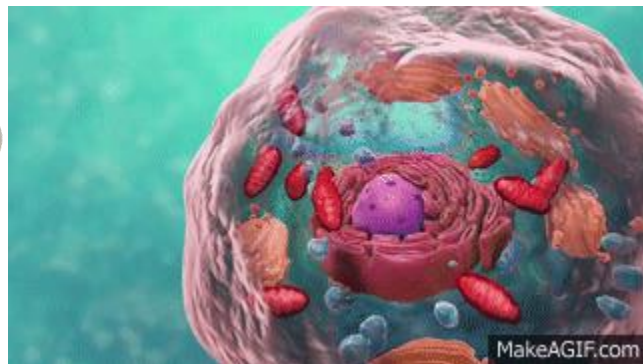
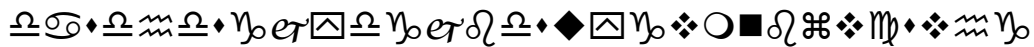


a biotechne® brand

ACD Document Types:

- a. Manual
- b. SDS
- c. CofA
- d. Datasheet
- e. Protocol
- f. Supplementary Data

Brand	Document Type	Document Name	File Type
RND	Manual	RND-Doc123	PDF
Novus Biological	CofA	NB%^_Doc-456	XML
ACD	SDS	7526 ACD_ &*!	DOC



<https://resources.novusbio.com/manual/Manual-NBP2-62174-47378557.pdf>