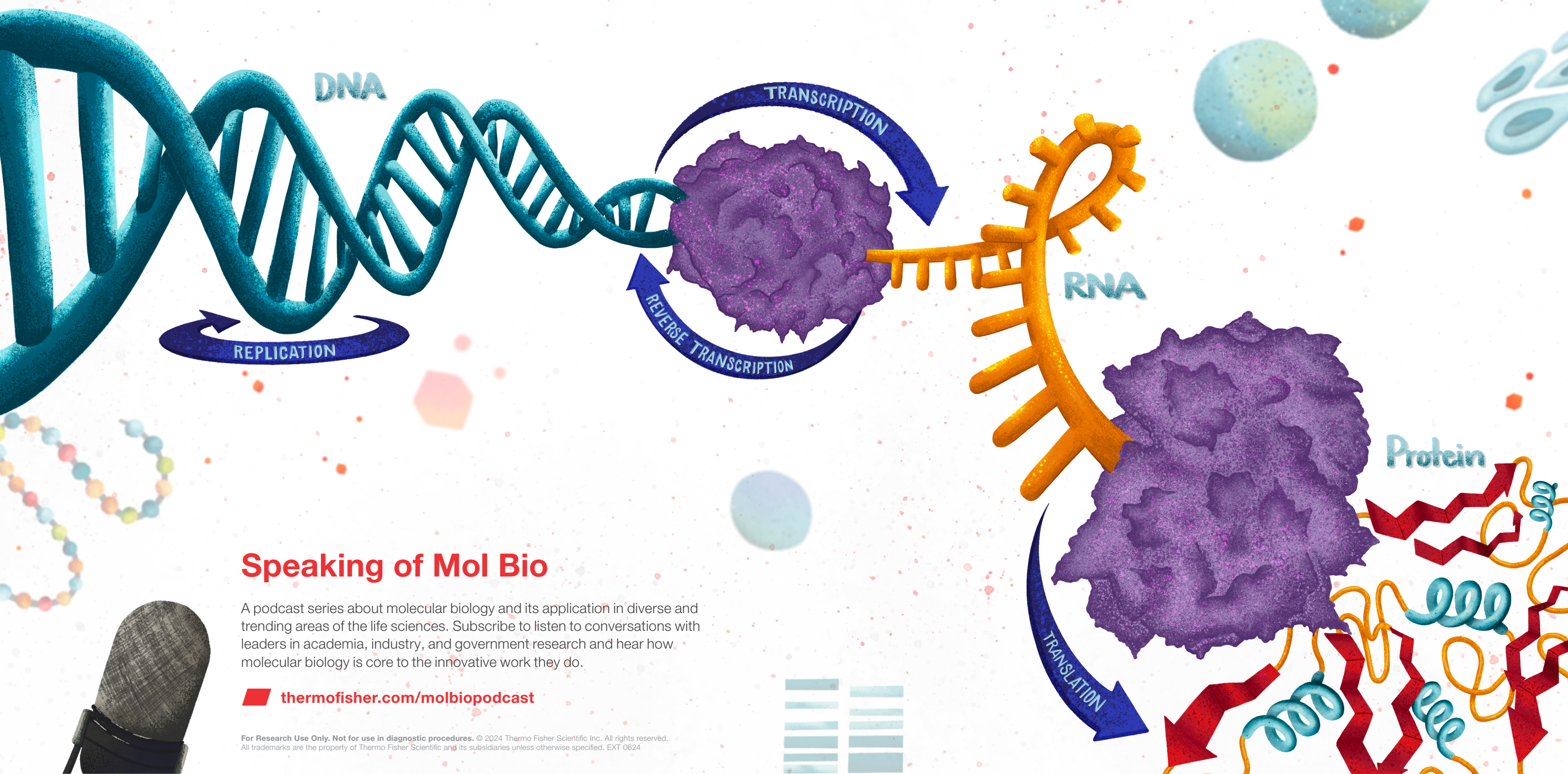


Molecular biology at the core



Speaking of Mol Bio

A podcast series about molecular biology and its application in diverse and trending areas of the life sciences. Subscribe to listen to conversations with leaders in academia, industry, and government research and hear how molecular biology is core to the innovative work they do.

 thermofisher.com/molbiopodcast