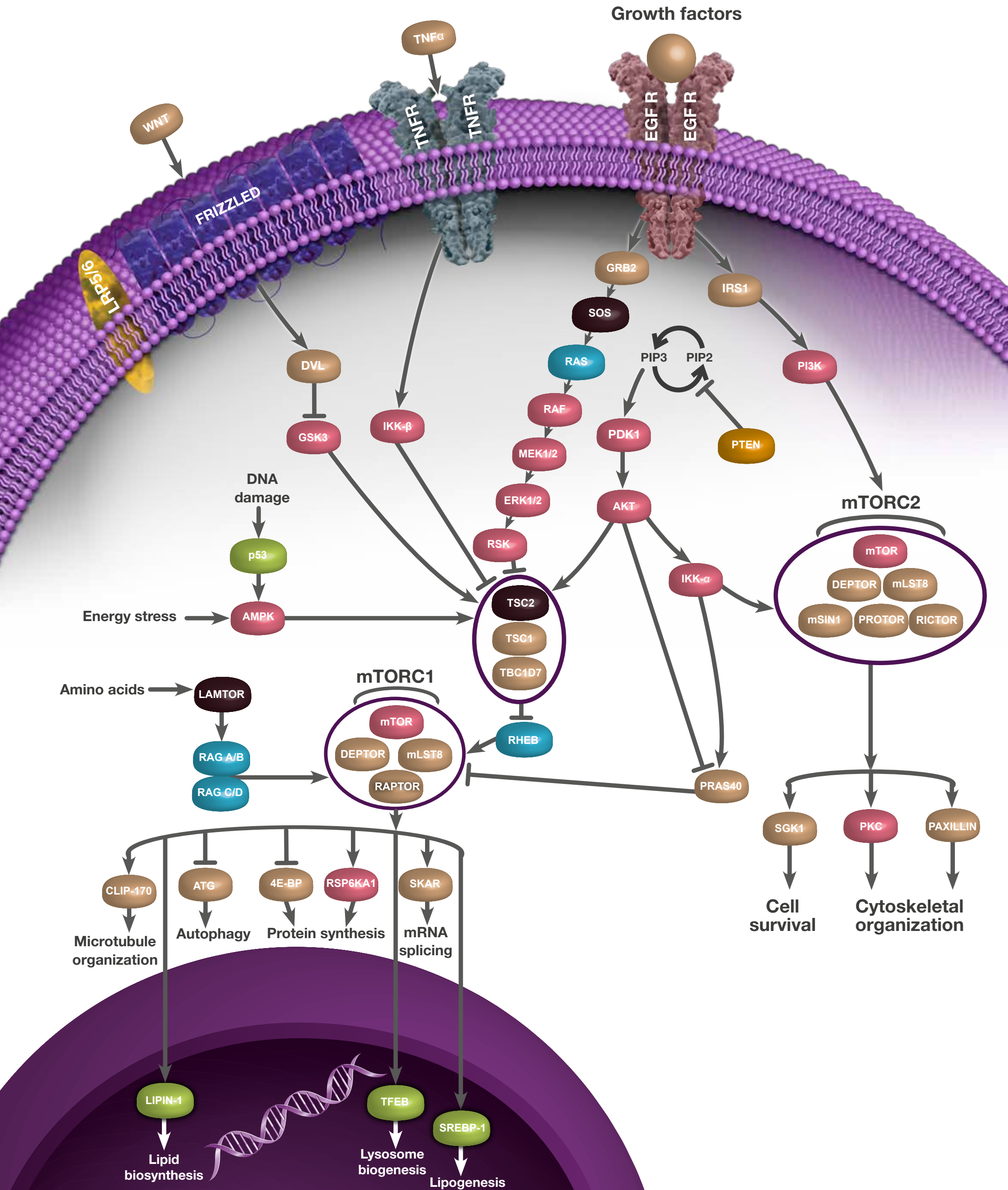


mTOR Signaling Pathway



BIO-RAD

Think Antibodies.
Think Bio-Rad.

