

PATHWAYS IN HUMAN CANCER

Revised Edition

The Pathways in Human Cancer guide summarizes some of the key signaling pathways involved in human cancer. It is the first guide to be published in human cancer, designed to be used by researchers in human cancer research. The guide provides a comprehensive overview of the pathways involved in human cancer, and is the first guide to be published in human cancer. The guide provides a comprehensive overview of the pathways involved in human cancer, and is the first guide to be published in human cancer.

The guide was developed by a team of experts in the field of human cancer research. It is the first guide to be published in human cancer, and is the first guide to be published in human cancer. The guide provides a comprehensive overview of the pathways involved in human cancer, and is the first guide to be published in human cancer.

<ul style="list-style-type: none"> Red box: protein-protein interaction Blue box: protein-DNA interaction Green box: protein-RNA interaction Yellow box: protein-lipid interaction Pink box: protein-carbohydrate interaction Grey box: protein-metal ion interaction 	<ul style="list-style-type: none"> Red arrow: activation Blue arrow: inhibition Green arrow: activation Yellow arrow: inhibition Pink arrow: activation Grey arrow: inhibition
---	--

