

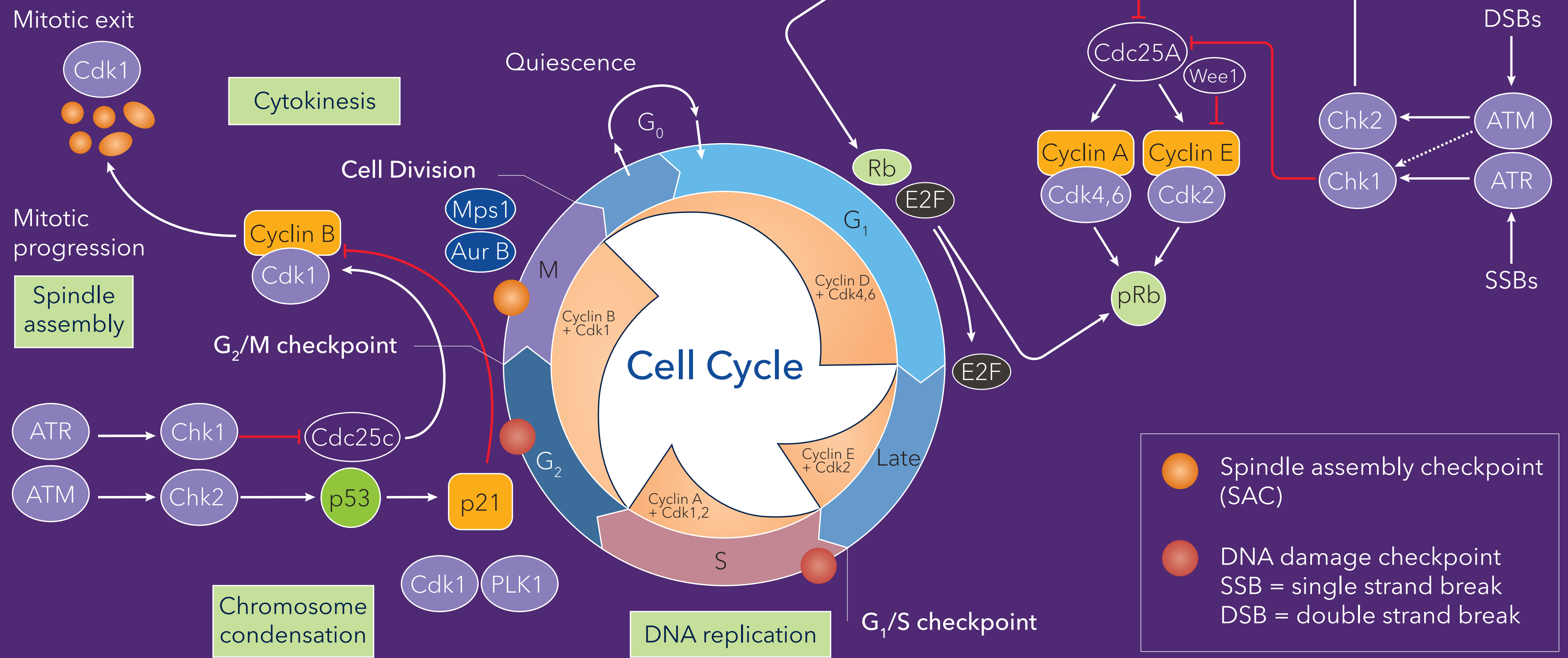
Cell Cycle and DNA Damage Repair

In normal cells, each stage of the cell cycle is tightly regulated. In cancer cells, many genes and proteins that influence the progression of the cell cycle are mutated or overexpressed - they become oncogenes. The proteins/molecules involved in the regulation of the cell cycle, in particular those with a role in DNA replication and DNA damage, are important cancer therapeutic targets.

Cell Cycle Progression and DNA Repair

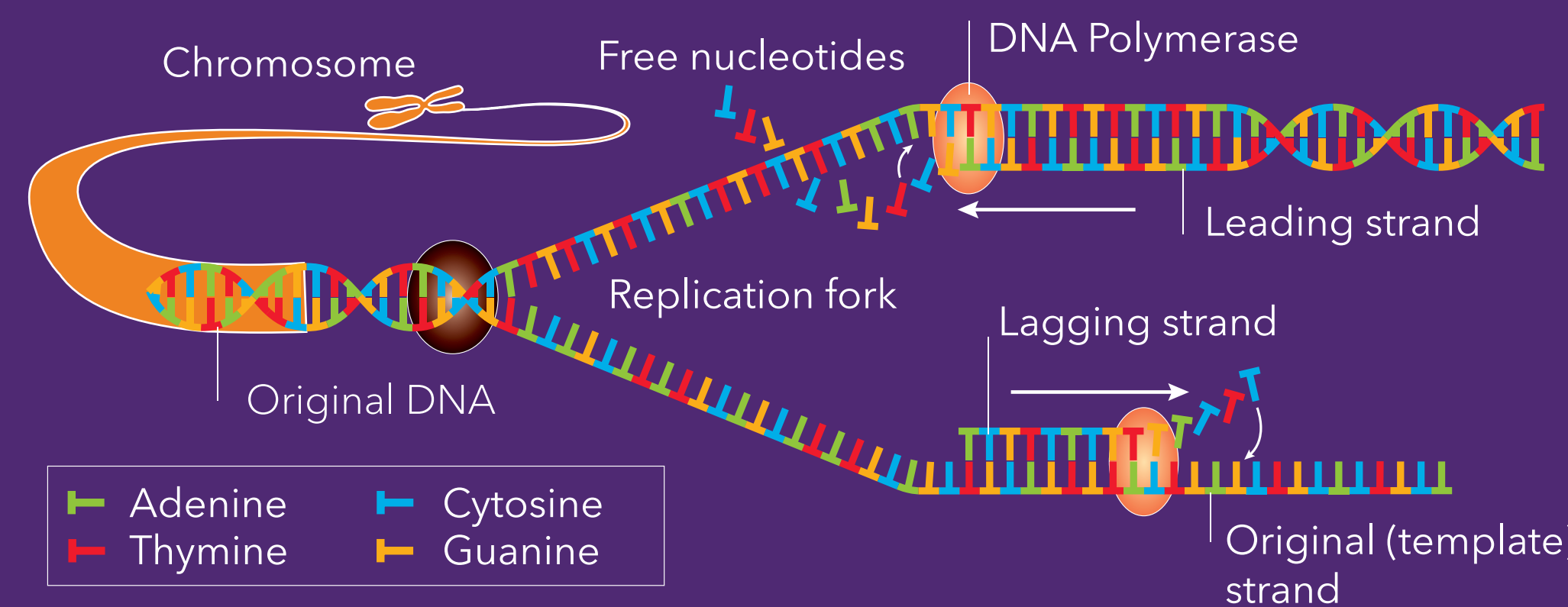
There are three major regulatory cell cycle checkpoints - G₁/S, intra-S phase and G₂/M phase. A cell can only pass through these checkpoints in the presence of stimulatory signals and in the absence of DNA damage. The cell cycle checkpoints are controlled by tumor suppressors and cyclin-dependent kinases (cdks). Cdks act in concert with their regulatory subunits, cyclins, to control cell cycle progression. Cdks are constitutively expressed and are regulated by several kinases and phosphatases, including Wee1 kinase and Cdc25 phosphatase.

At specific points in the cell cycle, DNA damage is detected and repaired. This process is initiated by the DNA damage sensors, ATM and ATR kinase. Checkpoint kinases Chk1 and Chk2 initiate signaling cascades that activate DNA damage checkpoints in G₁ and G₂. The spindle assembly checkpoint (SAC) delays anaphase of mitosis until all chromosomes are properly aligned on the spindle, preventing aneuploidy. Kinases including aurora kinase B (Aur B), PLK1 and Mps1 are implicated at various control points in the cell cycle.



DNA Replication

DNA replication occurs in five stages during S-phase; initiation, unwinding, primer synthesis, elongation and termination. Helicase enzymes "unwind" the DNA double helix, and telomerases reduce the resulting torsional strain, the single strands are now exposed and the replication fork is initiated. The leading strand of DNA is synthesized by Pol ε and the lagging strand is synthesized by Pol δ. PCNA is a cofactor for both DNA polymerase δ and ε, where it acts as a DNA clamp, which is important in both DNA synthesis and repair. At the end of the termination phase, DNA ligases form a phosphodiester bond, which joins the DNA strands together, forming new doubled stranded DNA.



Tocris Products

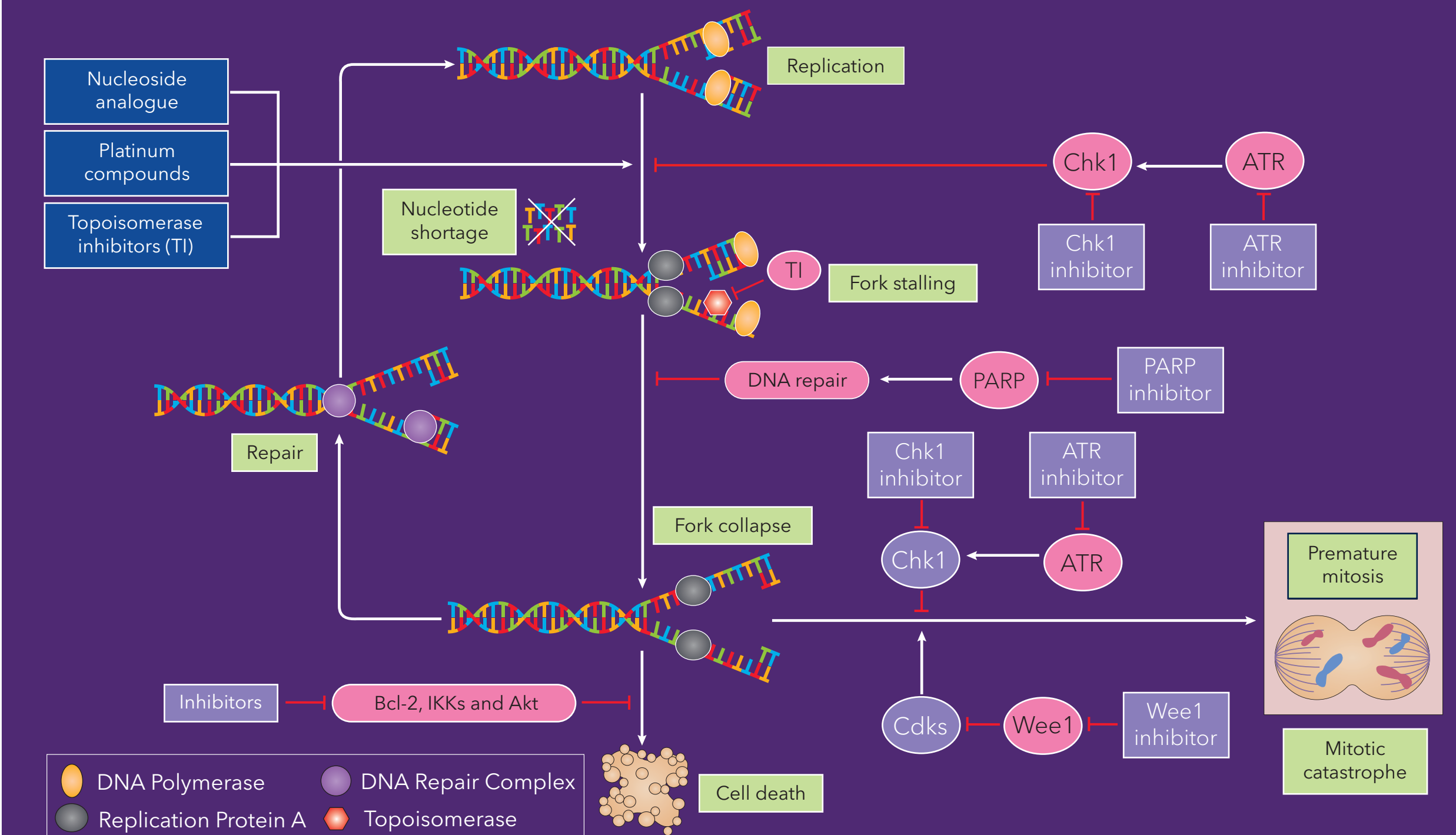
ATM & ATR Kinase AZ 20 AZ 5704 KU 55933 Mirin	Casein Kinase 2 TBB TBCA TTP 22	Cyclin-dependent Kinase BSJ-03-123 BSJ-04-132 CDK8/19i FMF-04-159-2 Kenpaullone NVP 2 PD 0332991 Purvalanol A Purvalanol B Ro 3306 Senexin A THAL SNS 032	Hsp70 VER 155008	p53 Nutlin-3 PRIMA-1MET RITA
Aurora Kinase Hesperadin TC-A 2317 ZM 447439	Cdc25 Phosphatase NSC 663284 NSC 95397	Cell Cycle Inhibitors 10058-F4 Methotrexate Narciclasine Pyridostatin	Hsp90 17-AAG	Polo-like Kinase GW 843682X TC-S
Calpain Acetyl-Calpastatin (184-210) (human) E 64 MG 132 PD 150606	Checkpoint Kinase CCT 241533 LY 2603618 NSC 109555 PD 407824 PF 477736 SB 218078 TCS 2312	DNA-dependent Protein Kinase NU 7026 NU 7441	Kinesin Dimethylenastron K 858 Monastrol S-Trityl-L-cysteine	Poly (ADP-ribose) Polymerase Nicotinamide PJ 34 WIKI 4 XAV 939
Casein Kinase 1 D 4476 PF 4800567 PF 5006739 PF 670462	DNA, RNA and Protein Synthesis L189 Mithramycin A NSC 617145	DNA, RNA and Protein Synthesis L189 Mithramycin A NSC 617145	Microtubules Docetaxel Flutax 1 Taxol Vinblastine Vincristine	Telomerase BRACO 19 BIBR 1532 RHPS 4 TMPyP4 tosylate

References

Annunziata and O'Shaughnessy (2010) *Clin. Cancer Res.* 16 4517.
 Barr et al (2004) *Nat. Rev. Mol. Cell Biol.* 5 429.
 Dobbstein and Sorensen (2015) *Nat. Rev. Drug Discov.* 14 405.
 Fu et al (2012) *Nat. Rev. Cancer* 12 104.
 Hochegger et al (2008) *Nat. Rev. Mol. Cell Biol.* 9 910.
 Lapenna and Giordano (2009) *Nat. Rev. Drug Discov.* 8 547.
 Lord and Ashworth (2012) *Nature* 481 287.
 Malumbres and Barbacid (2009) *Nat. Rev. Cancer.* 9 153.
 Rastogi and Mishra (2012) *Cell Div.* 7 26.
 Williams and Stoeber (2012) *Pathol.* 226 352.

Targeting Cancer Cells

Enhancing replicative stress by targeting critical DNA replication checkpoints and replication machinery, as well as depleting nucleotides, encourages fork stalling and fork collapse, which leads to mitotic catastrophe and death in cancer cells.



NOTE: This poster conveys a general overview and should be considered neither comprehensive nor definitive. The details of this information are understood to be subject to interpretation.

Learn more | tocris.com