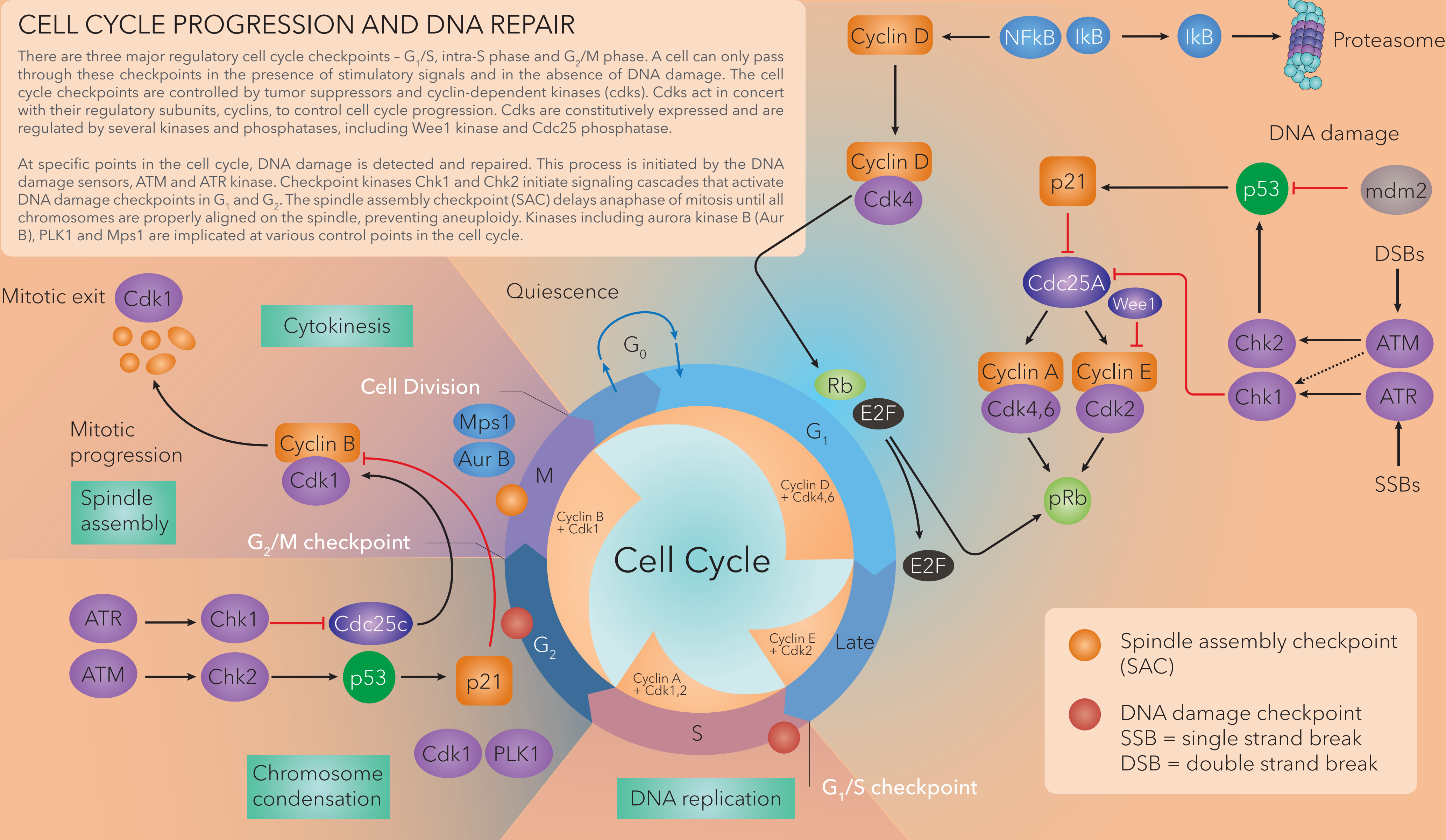


In normal cells, each stage of the cell cycle is tightly regulated. In cancer cells, many genes and proteins that influence the progression of the cell cycle are mutated or overexpressed – they become oncogenes. The proteins/ molecules involved in the regulation of the cell cycle, in particular those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



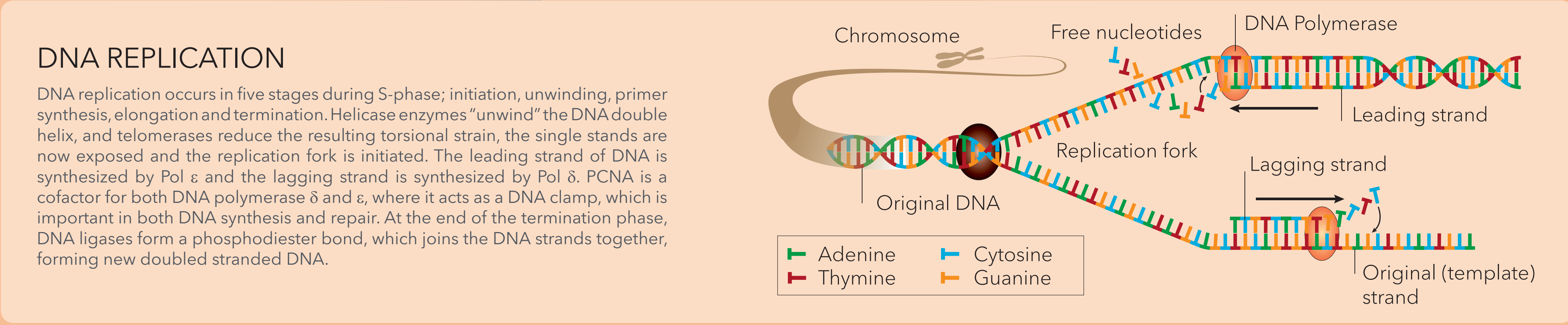
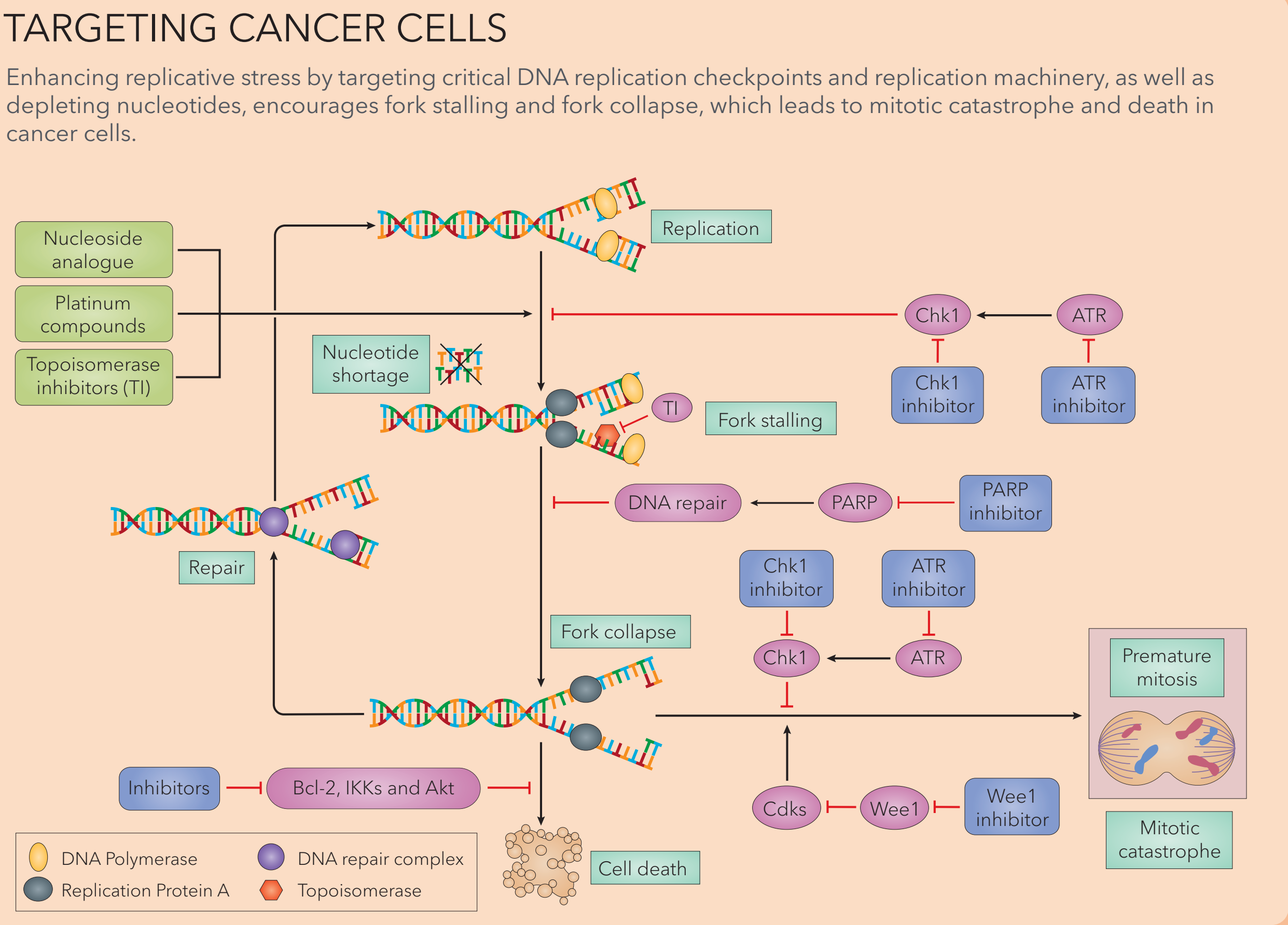
PRODUCTS AVAILABLE FROM TOCRIS

ATM & ATR Kinase	AZ 20, AZ 5704, KU 55933, Mirin, VE 821	DNA, RNA and Protein Synthesis	4E1RCat, L189, Mithramycin A, NSC 617145
Aurora Kinases	Hesperadin, TC-A 2317, VX 680, ZM 447439	Hsp70	VER 155008
Calpains	Acetyl-Calpastatin (184-210) (human), Calpeptin, E 64, MG 132, PD 150606	Hsp90	17-AAG
Casein Kinase 1	D 4476, IC 261, PF 4800567, PF 5006739, PF 670462	IRE1	APY 29
Casein Kinase 2	TBB, TBCA, TMCB, TTP 22	Kinesin	Dimethylnastron, K 858, Monastrol, SB 743921, S-Trityl-L-cysteine
Cdc25 Phosphatase	NSC 663284, NSC 95397	Microtubules	Docetaxel, Dolastatin 10, Flutax 1, Taxol, Vinblastine, Vincristine
Cell Cycle Inhibitors	10058-F4, Methotrexate, Narciclasine, Pyridostatin pentahydrochloride	Monopolar Spindle 1 Kinase	AZ 3146
Checkpoint Kinases	CCT 241533, LY 2603618, NSC 109555, PD 407824, PF 477736, SB 218078, TCS 2312	p53	Nutlin-3, PRIMA-1MET, RITA
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	Polo-like Kinase	GW 843682X, TC-S 7005
DNA-dependent Protein Kinase	NU 7026, NU 7441	Poly (ADP-ribose) Polymerase	Nicotinamide, PJ 34, WIKI 4, XAV 939
		Telomerase	BRACO 19 trihydrochloride, BIBR 1532, RHPS methosulfate, TMPyP4 tosylate

REFERENCES

Annunziata and O'Shaughnessy (2010) *Clin. Cancer Res.* 16 4517.
Barr et al (2004) *Nat. Rev. Mol. Cell Biol.* 5 429.
Dobbelstein 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 Stoerber (2012) *Pathol.* 226 352.



To request a copy of the Tocris Cancer Product Guide, or to view the PDF, please visit | [toocris.com](https://www.tocris.com)

Adapted from Edition 3 of the Tocris Cancer Product Guide