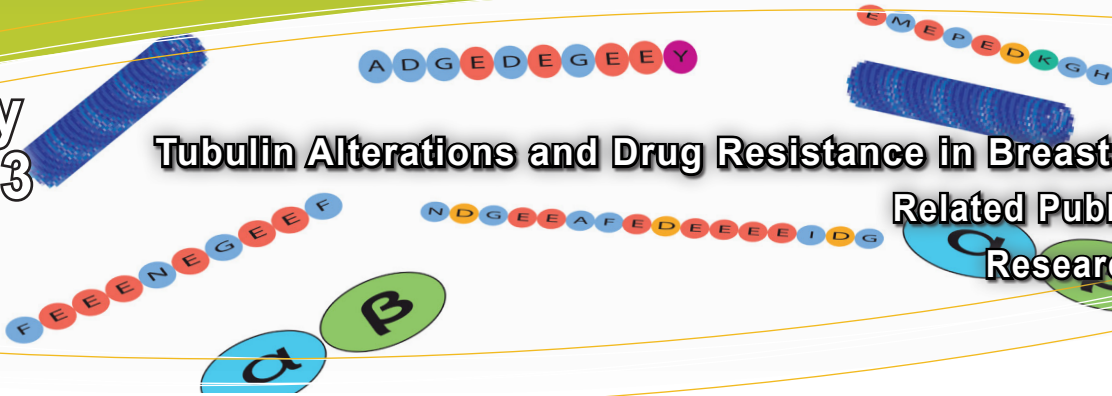




Helping advance science,  
one protein at a time.

May  
2023

## Tubulin Alterations and Drug Resistance in Breast Cancer Related Publications Research Tools



### Sponsored Conferences

JCS Imaging Cell  
Dynamics  
Pestana Palace Hotel,  
Lisbon Portugal  
May 14<sup>th</sup>-17<sup>th</sup>

Gordon Scientific  
Phagocytes  
Waterville Valley, NH  
June 3<sup>rd</sup> - 4<sup>th</sup>

### Cytoskeleton Products

- Actin Proteins
- Activation Assays
- Antibodies
- ECM Proteins
- ELISA Kits
- G-LISA® Kits
- Live Cell Imaging
- Pull-down Assays
- Motor Proteins
- Small G-Proteins
- Tubulin & FtsZ Proteins

### Contact Us

P: 1 (303) 322.2254  
F: 1 (303) 322.2257  
E: [cserve@cytoskeleton.com](mailto:cserve@cytoskeleton.com)  
W: [cytoskeleton.com](http://cytoskeleton.com)

### For More News

W: [cytoskeleton.com /blog/](http://cytoskeleton.com/blog/)

## Tubulin Alterations and Drug Resistance in Breast Cancer

### Introduction

Breast cancer (BC) is currently accountable for 1 in 8 cancer diagnoses worldwide, with 2.3 million new patients annually, both sexes combined.<sup>1</sup> In women, it constitutes a quarter of all cancer cases and has become the most frequently detected form of the disease in 2020. This matter has been escalating globally, especially in developing countries.<sup>2</sup> In the United States, mortality decreased 41% since 1989; however, the descending trend has slowed, stressing the need for enhanced treatments.<sup>3,4</sup> During diagnosis, BC is classified according to the expression of hormone receptors (estrogen and progesterone receptors, primarily), the overexpression of the human epidermal growth factor receptor 2 (oncogene HER2),<sup>5</sup> and with a pathological classification that evaluates nuclear pleomorphism, mitotic count, and tubule count. Additionally, diverse alterations in the microtubule network have been detected and characterized in BC, including irregular expression of tubulin isotypes and abnormal tubulin post-translational modifications.<sup>6</sup> Microtubules are protein structures that constitute dynamic pillars of the cytoskeleton and have major roles in key cellular processes, such as mitotic chromosome segregation, cell shape and motility, and inner cellular transport.<sup>7</sup> Alterations in these essential structures in BC are associated with poor prognosis and response to treatment, and a more aggressive disease.<sup>6</sup>

### Molecular Markers Used for Drug Response Assessment

Molecular markers expressed by cancerous cells are mutated/modified proteins that bind to hormones, gene expression patterns, and altered DNA sequences that can function as indicators of response to specific therapies.<sup>8</sup> The underlying cause of this phenomenon relies on DNA mutations that vary from person to person and can affect drug efficacy. Among the most well-known molecular markers related to drug resistance, there is the oncogene HER2, whose expression is amplified in 20-30% of BC cases and is regarded as a marker of poor prognosis. HER2 overexpression is linked to resistance to antihormonal and cytotoxic therapies.<sup>8</sup> A further example is the low transcriptional expression of CYP2D6, which predicts resistance to chemotherapy with tamoxifen (a drug that blocks estrogen stimulation) in BC.<sup>8,9</sup> In some cases, drug resistance has hampered the clinical success of microtubule-targeting agents (MTAs) in BC (and other cancers). More recent evidence has indicated that specific tubulin isotypes, such as class III  $\beta$ -tubulin, can leave microtubule drugs without effect in cells. As a result, tubulin isotypes are currently under study as potential prognostic biomarkers.<sup>10</sup>

### Historical Use of Microtubule-Targeting Agents (MTAs) for Cancer Therapy

Given the central role of microtubules in vital cellular processes, agents that target these structures can impair normal cell function and will often lead to cell death. As cancer cells divide rapidly, they are more susceptible to cell cycle arrest-induced death; thus, MTAs have been counted among the cancer treatments of choice for decades. However, MTAs with better-targeting abilities are still sought after. The first MTA to be introduced in clinical cancer therapy was paclitaxel, which has been used for BC since 1994. Subsequent MTAs that were approved for BC treatment were the semi-synthetic taxane docetaxel, the more recent taxanes larotaxel and ixabepilone, and the vinca alkaloids.<sup>11-14</sup> In the present day, paclitaxel and vinca alkaloids have been established as the standard drugs in the management of different cancers, including BC.<sup>14</sup> Nonetheless, drug resistance due to long-term use and solubility problems have incentivized the development of novel MTAs for BC treatment, including epothilone, eribulin, auristatin, and maytansine.<sup>15-17</sup> The consequences of MTA therapeutic use on microtubule dynamics entail disruption of intracellular cell transport, cessation of cell division, and triggering of cell death. Noteworthy, the vast majority of MTAs interfere with microtubule functions by acting on their  $\beta$ -tubulin subunit.<sup>18</sup>

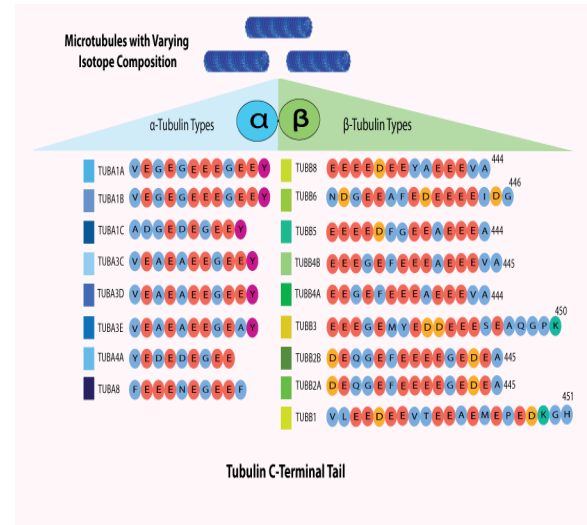


Figure 1. Schematic of alpha and beta tubulin isotypes. Shown are the varying C-terminal sequences of these isotypes.



Helping advance science,  
one protein at a time.

# Tubulin PRODUCTS

## Continued from Page 1

### Recent Findings: Alterations in $\beta$ III-tubulin Gene Expression are Associated with MTA Resistance in BC Patients

In mammals, cells are known to express not less than eight different  $\beta$ -tubulin isotypes, identified as  $\beta$ I,  $\beta$ IIa,  $\beta$ IIb,  $\beta$ III,  $\beta$ IVa,  $\beta$ IVb,  $\beta$ V, and  $\beta$ VI (Figure 1).<sup>19-20</sup> Thorough research has evidenced that alterations in some of these isotypes in cancer cells were associated with resistance to MTAs.<sup>10,21,22</sup> In particular, an increased abundance of  $\beta$ III-tubulin represents the most prevalent mechanism concerning MTAs resistance in various tumor types, including BC.<sup>10,21</sup> In this regard, Lopus et al evaluated the effects of MTA ixabepilone on BC cells with and without the *in vitro* removal and knockdown of  $\beta$ III-tubulin.<sup>23</sup> They concluded that  $\beta$ III-tubulin expression inhibits the antitumor effects of ixabepilone, denoting that increased  $\beta$ III-tubulin could contribute to ixabepilone resistance. Scherbakov et al analyzed the impact of long-term incubation of ER $\alpha$ -positive BC cells with docetaxel.<sup>24</sup> For this purpose, they evaluated the expression of signaling proteins by immunoblotting and flow cytometry and assessed ER $\alpha$  activity via gene reporter assay. They found that the cells with the highest  $\beta$ III-tubulin levels were resistant to docetaxel, while those with the lowest  $\beta$ III-tubulin expression were sensitive to the drug. Extensive research in this field continues today, and further studies have claimed that  $\beta$ III-tubulin does not work on its own.<sup>25</sup> Moreover, it has been found that the molecular pathways influenced by  $\beta$ III-tubulin depend on the cell and cancer type, which could explain why some tumor types do not show poor results and resistance to MTAs at high  $\beta$ III-tubulin levels.<sup>25</sup>

### Summary and Future Perspectives

In a global context of high rates of BC, it is of utmost relevance to explore, detect, and study genetic biomarkers that can predict individual responses to the administered drugs. Overexpression of  $\beta$ III-tubulin has been found in numerous cancer types, including BC, and it has been linked to poor response and resistance to various MTA. As a result,  $\beta$ III-tubulin has become a considerable biomarker candidate. Nevertheless, for this knowledge to be fully exploited in clinical medicine, further conclusive research is needed, together with more studies considering the particular contexts of the patients and rigorous clinical evaluation.

## Purified Tubulin

Product	Amount	Cat #
<b>Tubulin protein (&gt;99% pure)</b> Source: porcine brain	1 x 1 mg 5 x 1 mg	T240-A T240-B

## Tubulin Proteins

Product	Amount	Cat #
<b>Microtubule associated protein rich fraction</b> Source: porcine brain	1 x 100 $\mu$ g 5 x 100 $\mu$ g	MAPF-A MAPF-C
<b>Tubulin protein (&gt;99% pure)</b> Source: porcine brain	1 x 1 mg 5 x 1 mg	T240-A T240-B
<b>Acetylated Tubulin Protein Source: Porcine Brain</b> Source: porcine brain	1 x 500 $\mu$ g	TAC01

## Tubulin Live Cell Imaging Products

Product	Ex / Em	Amount	Cat #
<b>SiR-Tubulin™ Kit</b> Includes SiR-Tubulin, and Verapamil	630 / 680 nm	50 nmol	CY-SC002
<b>SiR700-Tubulin Kit</b> 35 nmol SiR700-Tubulin and 1 $\mu$ mol verapamil	680 / 720 nm	50 nmol	CY-SC014
<b>SPY555-Tubulin Kit</b> Includes SPY555-Tubulin and Verapamil	555 / 580 nm	100 stains	CY-SC203
<b>SPY650-Tubulin</b> Includes SPY650-Tubulin and Verapamil	652 / 674 nm	100 stains	CY-SC503

## References

- Arnold M, Morgan E, Rungay H, Mafra A, Singh D, Laversanne M, et al. Current and future burden of breast cancer: Global statistics for 2020 and 2040. *The Breast*. 2022 Dec 1;66:15–23.
- Heer E, Harper A, Escandor N, Sung H, McCormack V, Fidler-Benaoudia MM. Global burden and trends in premenopausal and postmenopausal breast cancer: a population-based study. *Lancet Glob Heal*. 2020 Aug 1;8(8):e1027–37.
- Siegel RL, Miller KD, Fuchs HE, Jemal A. *Cancer Statistics, 2021*. *CA Cancer J Clin* [Internet]. 2021 Jan [cited 2023 Mar 15];71(1):7–33. Available from: <https://pubmed.ncbi.nlm.nih.gov/33433946/>
- American Cancer Society. *Breast Cancer Facts & Figures 2019-2020*. Atlanta; 2019.
- Li C, Fan Z, Lin X, Cao M, Song F, Song F. Parity and risk of developing breast cancer according to tumor subtype: A systematic review and meta-analysis. *Cancer Epidemiol*. 2021 Dec 1;75:102050.
- Parker AL, Kavallaris M, McCarroll JA. Microtubules and Their Role in Cellular Stress in Cancer. *Front Oncol* [Internet]. 2014 [cited 2023 Apr 4];4. Available from: <https://pmc/articles/PMC4061531/>
- Logan CM, Menko AS. Microtubules: Evolving roles and critical cellular interactions. *Exp Biol Med* (Maywood) [Internet]. 2019 Nov 1 [cited 2023 Mar 20];244(15):1240–54. Available from: <https://pubmed.ncbi.nlm.nih.gov/31387376/>
- Banin Hirata BK, Oda JMM, Losi Guembarovski R, Ariza CB, Oliveira CEC De, Watanabe MAE. Molecular markers for breast cancer: prediction on tumor behavior. *Dis Markers* [Internet]. 2014 [cited 2023 Mar 22];2014. Available from: <https://pubmed.ncbi.nlm.nih.gov/24591761/>
- Mehta S, Lasham A, Blenkiron C, Shelling A, Muthukaruppan A, Laking G, et al. Predictive and prognostic molecular markers for cancer medicine. *Ther Adv Med Oncol* [Internet]. 2010 [cited 2023 Mar 22];2(2):125. Available from: <https://pmc/articles/PMC3126011/>
- Parker AL, Teo WS, McCarroll JA, Kavallaris M. An Emerging Role for Tubulin Isotypes in Modulating Cancer Biology and Chemotherapy Resistance. *Int J Mol Sci* [Internet]. 2017 Jul 4 [cited 2023 Mar 24];18(7). Available from: <https://pubmed.ncbi.nlm.nih.gov/28677634/>
- Engels FK, Sparreboom A, Mathot RAA, Verweij J. Potential for improvement of docetaxel-based chemotherapy: a pharmacological review. *Br J Cancer* [Internet]. 2005 Jul 7 [cited 2023 Apr 5];93(2):173. Available from: <https://pmc/articles/PMC2361544/>
- Ren S, Wang Y, Wang J, Gao D, Zhang M, Ding N, et al. Synthesis and biological evaluation of novel larotaxel analogues. *Eur J Med Chem*. 2018 Aug 5;156:692–710.
- Rugo HS, Barry WT, Moreno-Aspitia A, Lyss AP, Cirrincione C, Leung E, et al. Randomized Phase III Trial of Paclitaxel Once Per Week Compared With Nanoparticle Albumin-Bound Nab-Paclitaxel Once Per Week or Ixabepilone With Bevacizumab As First-Line Chemotherapy for Locally Recurrent or Metastatic Breast Cancer: CALGB 40502/NCCTG N063H (Alliance). *J Clin Oncol* [Internet]. 2015 Jul 7 [cited 2023 Apr 5];33(21):2361. Available from: <https://pmc/articles/PMC4500830/>
- Cermák V, Dostál V, Jelínek M, Libusová L, Kovář J, Rösel D, et al. Microtubule-targeting agents and their impact on cancer treatment. *Eur J Cell Biol* [Internet]. 2020 May 1 [cited 2023 Mar 20];99(4). Available from: <https://pubmed.ncbi.nlm.nih.gov/32414588/>
- Lambert JM, Charl RVJ. Ado-trastuzumab Emtrastine (T-DM1): an antibody-drug conjugate (ADC) for HER2-positive breast cancer. *J Med Chem* [Internet]. 2014 [cited 2023 Apr 5];57(16):6949–64. Available from: <https://pubmed.ncbi.nlm.nih.gov/24967516/>
- Perez-Garcia JM, Cortes J. The safety of eribulin for the treatment of metastatic breast cancer. <https://doi.org/10.1080/1474033820191608946> [Internet]. 2019 May 4 [cited 2023 Apr 5];18(5):347–55. Available from: <https://www.tandfonline.com/doi/abs/10.1080/14740338.2019.1608946>
- Chung SW, Cho YS, Choi JU, Kim HR, Won TH, Kim SY, et al. Highly potent monomethyl auristatin E prodrug activated by caspase-3 for the chemoradiotherapy of triple-negative breast cancer. *Biomaterials*. 2019 Feb 1;192:109–17.
- Steinmetz MO, Prota AE. Microtubule-Targeting Agents: Strategies To Hijack the Cytoskeleton. *Trends Cell Biol*. 2018 Oct 1;28(10):776–92.
- Gadadhar S, Bodakuntla S, Natarajan K, Janke C. The tubulin code at a glance. *J Cell Sci* [Internet]. 2017 Apr 15 [cited 2023 Mar 24];130(8):1347–53. Available from: <https://journals.biologists.com/jcs/article/130/8/1347/56730/The-tubulin-code-at-a-glance>
- Ludueña RF. A hypothesis on the origin and evolution of tubulin. *Int Rev Cell Mol Biol* [Internet]. 2013 [cited 2023 Mar 24];302:41–185. Available from: <https://pubmed.ncbi.nlm.nih.gov/23351710/>
- Kanakkanthara A, H. Teesdale-Spittle P, H. Miller J. Cytoskeletal Alterations that Confer Resistance to Anti-tubulin Chemotherapeutics. *Anticancer Agents Med Chem*. 2013 Feb 15;13(1):147–58.
- Prassanawar SS, Panda D. Tubulin heterogeneity regulates functions and dynamics of microtubules and plays a role in the development of drug resistance in cancer. *Biochem J* [Internet]. 2019 May 12 [cited 2023 Mar 24];476(9):1359–76. Available from: <https://pubmed.ncbi.nlm.nih.gov/31085712/>
- Lopus M, Smiyun G, Miller H, Oroudjev E, Wilson L, Jordan MA. Mechanism of action of ixabepilone and its interactions with the  $\beta$ III-tubulin isotype. *Cancer Chemother Pharmacol* [Internet]. 2015 Nov 1 [cited 2023 Apr 12];76(5):1013–24. Available from: <https://pubmed.ncbi.nlm.nih.gov/26416565/>
- Scherbakov AM, Basharina AA, Sorokin D V, Mikhaevich EI, Mizaeva IE, Mikhaylova AL, et al. Targeting hormone-resistant breast cancer cells with docetaxel: a look inside the resistance. *Cancer Drug Resist* [Internet]. 2023 Feb 7 [cited 2023 Mar 25];6(1):103–15. Available from: <https://cdrjournal.com/article/view/5423>
- Kanakkanthara A, Miller JH.  $\beta$ III-tubulin overexpression in cancer: Causes, consequences, and potential therapies. *Biochim Biophys Acta Rev Cancer* [Internet]. 2021 Dec 1 [cited 2023 Mar 25];1876(2). Available from: <https://pubmed.ncbi.nlm.nih.gov/34364992/>

## Tubulin Kits

Product	Assays	Cat #
<b>Tubulin polymerization HTS assay using &gt;97% pure tubulin</b> OD based - Porcine	24	BK004P
<b>Tubulin polymerization assay using &gt;99% pure tubulin</b> OD based - Porcine	24-30	BK006P
<b>Tubulin polymerization assay using &gt;99% pure tubulin</b> Fluorescence based	96	BK011P
<b>Microtubule Binding Protein Spin-Down Assay Biochem Kit</b>	50-100	BK029
<b>Microtubule/Tubulin In Vivo Assay Biochem Kit</b>	30-100	BK038